



Green Fluorescent GECO Calcium (Ca²⁺) Assay

July 6, 2026

Product Info & Protocol

About these Assays (Table 1)	2
Assay Kit Materials and Storage (Table 2)	3
Additional Materials Required	3
Biosafety Considerations	3
Terms of Sale	3
Protocol for Use	4
Before your First Experiment – Optimize by Titrating the Sensor	4
Day 1 – Transduce and Plate Cells	5
Suggestions for Assays in Adherent Cells	6
Scaling for 384 Wells	6
Alternative Spinoculation Protocol	7
Day 2 – Measuring Ca ²⁺ Response	8
Measure Fluorescence on a Plate Reader or Imaging System	8
Fluorescence Detection (Tables 3-5)	9
Fluorescence Properties (Fig. 1)	10
Timing (Fig. 2)	10
Data Analysis	10
Assay Optimization	11
Optimizing Fluorescence	11
Optimizing Expression for Your Cell Type	11
Use the Positive Control	11
Optimizing Receptor Expression	11
Troubleshooting Guide	12
GECO Ca ²⁺ Sensors – in the Literature	16
References	16
Related Products	16
Contact Us	16



About these Assays

The green fluorescent Ca²⁺ sensor, Green GECO, increases in fluorescence intensity in response to increases in intracellular calcium. Green GECO was developed in [Robert Campbell's lab at the University of Alberta \[1\]](#). Changes in fluorescence in living cells can be detected with epifluorescence microscopy or on automated fluorescence plate readers. This green GECO calcium sensor can be combined with different colored sensors to measure multiple signals simultaneously. For example, product #U0650G, Green GECO Ca²⁺ sensor can be combined with a red fluorescent sensor for DAG (product # D0350R) to better understand the kinetics of coordinated, parallel signaling processes in living cells [5].

For sensitive and/or difficult to transduce cell types, such as primary neurons, we offer purified, high-titer BacMam stock. Please [contact us](#) to discuss whether purified BacMam is the right product for you.

The GECO Ca²⁺ Assay protocol is optimized for measuring Ca²⁺ responses in rapidly dividing, immortalized cell lines on a 96-well plate, and has been validated in live HEK 293T cells [3]. Green GECO users have [published results](#) in a wide variety of cell types. This assay is very robust and can be used for live-cell imaging or for screening on automated fluorescence plate readers. For use in iPSC-derived or adherent cells, see [Suggestions for Assays in Adherent Cells](#) section. For use in difficult to transduce cell types, see [Alternative Spinoculation Protocol](#) section.

The protocol steps are simple and will be performed over two consecutive days. Please remember to [optimize the assay for your conditions](#) before performing your first experiment.

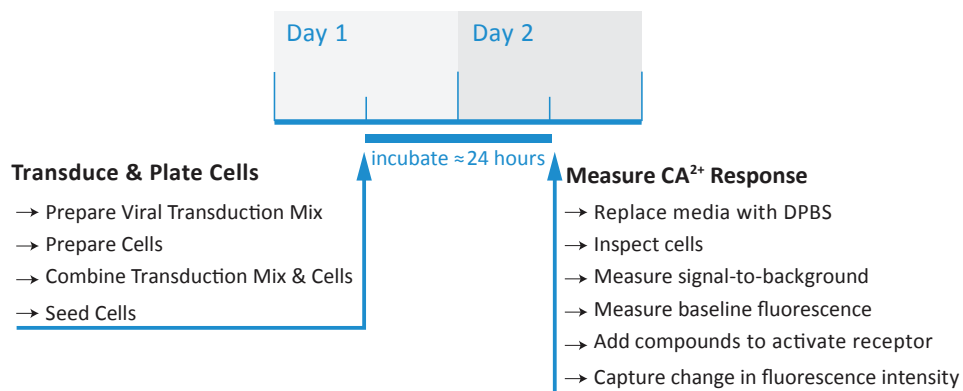
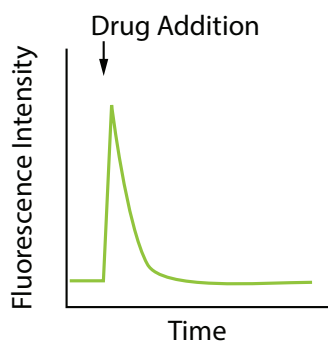
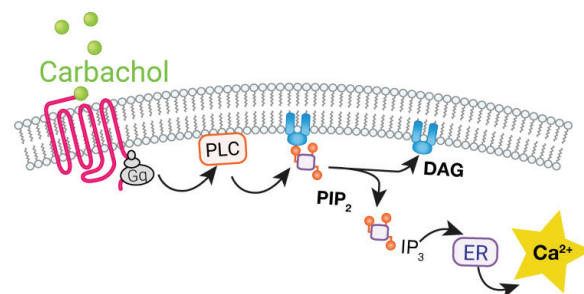


Table 1. Green GECO Calcium Kits

Product	Promoter	BacMam Version	Description	Recommended Use
#U0650G	CMV	Big Sky	Green GECO Ca ²⁺	Fluorescence imaging and plate reader assay
#U0600G	CMV	OG	Green Up G-GECO Ca ²⁺	Fluorescence imaging and plate reader assay (Z' < 0.5)
#U0608G	Ef1-α	OG	Green Up G-GECO Ca ²⁺	Fluorescence imaging and plate reader assay (Z' < 0.5)



Questions?

Call us, we can help!

+1 406-200-8321

info@montanamolecular.com

Our goal is to make your workflow easy and reproducible.






We'd love to hear about your research.



We also have a Troubleshooting Guide

Assay Kit Materials and Storage

BacMam stocks should be stored at 4°C protected from light in the original package. Store control agonist at -20°C. Avoid repeated freeze/thaw cycles. We recommend re-testing BacMam stock after storing for more than 12 months. If your BacMam stock has been purified, use it within 30 days for best results. Store sodium butyrate at 4°C.

Table 2. Materials in Kit	Details	Storage
Green GECO Ca ²⁺ sensor BacMam ≅ 2 × 10 ¹⁰ VG/mL in ESF 921 Insect Culture Medium (Expression Systems, product #96-001-01)	Green fluorescent sensor that changes in fluorescence intensity in response to increases in Ca ²⁺ . VG/mL is the number of viral genes per milliliter (see Biosafety Considerations section).	4°C
 sodium butyrate (Sigma Aldrich product #B5887) 500 mM in H ₂ O	Sodium butyrate is added to the culture to maintain BacMam expression. Other HDAC inhibitors may work as well.	4°C
 M1 muscarinic acetylcholine receptor BacMam in ESF 921 Insect Culture Medium (Expression Systems, product #96-001-01)	A Gq-coupled receptor provided as a positive control for the purpose of assay optimization. Contains a separate red fluorescent protein that is targeted to the nucleus.	4°C
 carbachol (Sigma Aldrich product number C4382) 25 mM in H ₂ O	Carbachol can be used to stimulate Gq signaling through the positive control, the M1 muscarinic acetylcholine receptor.	-20°C

Additional Materials Required (not included in kit)

1. Black, clear bottom microplate coated with a cell attachment factor. We recommend the following plates; 96-well [Greiner Bio-One \(#655946\)](#), 96-well Thermo Fisher Scientific ([#152037](#)), 384-well [Greiner Bio-One: \(#781946\)](#), 384-well [Corning: \(#354663\)](#).
2. Dulbecco's Phosphate Buffered Saline with Ca²⁺ and Mg²⁺ (PBS) available from VWR [\[4\]](#).
3. Cells and cell media. We recommend media with low autofluorescence such as EMEM, McCoy's 5A, and F12K culture media.

Biosafety Considerations

The BacMam vector carrying the fluorescent biosensor in these assays is a modified baculovirus, used for delivery to, and expression in, a wide variety of mammalian cell types including primary cultures.

BacMam is a modified baculovirus, *Autographa californica*, AcMNPV. The natural host of baculovirus is larvae of the order *Lepidoptera*. The BacMam vector in the kit is produced in Sf9 insect cells and is pseudotyped to infect mammalian cells. In mammalian cells, the baculovirus genome is silent and it cannot replicate to produce new virus. While it should be handled carefully, in a sterile environment, it is classified as a Biosafety Level 1 (BSL-1) reagent [\[7\]](#).

Other types of viruses are quantified in terms of plaque forming units (PFU) in cells from the natural host. Since BacMam is modified to produce expression in mammalian cells, we quantify the virus by measuring viral genes (VG) per milliliter (mL). Viral samples are prepared to release viral genomic DNA, then multiple dilutions of the preparation are run in qPCR using primers that are specific to the VSVG gene in the BacMam genome. Results are compared against a standard curve to generate an average titer for each viral stock. Check the label on the tube to find VG/mL for your stock.

This product is for research use only and is not for use or sale in human or animal diagnostic or therapeutic products.

Terms of Sale

All materials in this kit are provided without warranty, express or implied. User is responsible for making sure product use complies with applicable regulations. No right to resell products or any components of these products is conveyed. Use of materials is restricted to the intended purpose described in this protocol. Reverse engineering or modification of materials is not permitted. User agrees to accept these Terms of Sale before using materials.



Protocol for Use

This protocol is optimized for use in HEK 293T cells, however, it can be adjusted for use with practically any cell type. **Take the time to optimize the assay for your cell type and your particular conditions.**

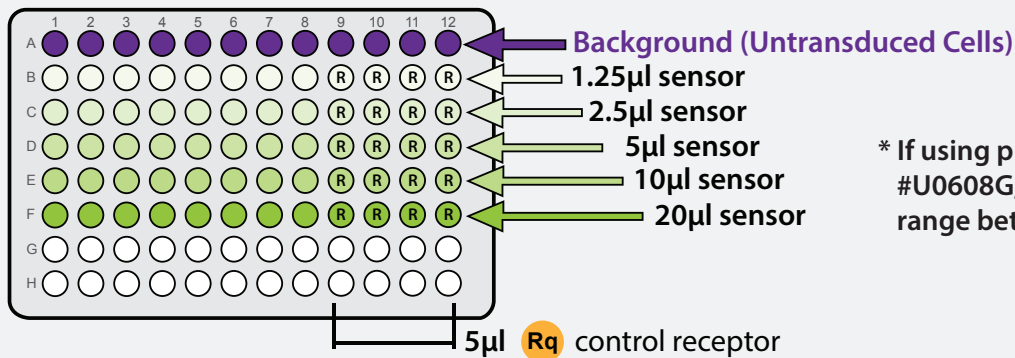
See our suggestions for [Assays in Adherent Cell Types](#), [Scaling for 384-well Plates](#), and [Alternative Spinoculation Protocol](#).

Before your first experiment: Optimize Your Assay by Titrating the Sensor

We recommend using a titration series to determine the best combination of signal above background, cell health, and sensor expression. Ideally, the signal in each well before drug addition should be at least 5 times above background.

Transduce and Plate Cells for Optimization

- Set up your plate. Be sure to include control wells (untransduced cells) in order to calculate signal-to-background. (For details on preparing cells and transduction mix, [refer to Day 1 procedure on page 5](#)).
- Perform titration to determine optimal sensor volume for your cells, as detailed in plate diagram and tables below.
- For each sensor volume, include a subset of wells that includes 5µl of Rq control receptor included in your kit. These wells will be used for your positive control.



* If using product #U0600G or #U0608G, we suggest a titration range between 2.5µL - 40µL.

Wells A1-A12 Background (with Untransduced Cells)

500 mM SB (sodium butyrate) 0.6µL

complete media 49.4µL

total volume per well 50µL

Wells B1-B8 Transduction Mix

GECO Ca²⁺ Sensor 1.25µL

500mM SB (sodium butyrate) 0.6µL

complete media 48.15µL

total volume per well 50µL

Wells B9-B12 Transduction Mix

GECO Ca²⁺ Sensor 1.25µL

500mM SB (sodium butyrate) 0.6µL

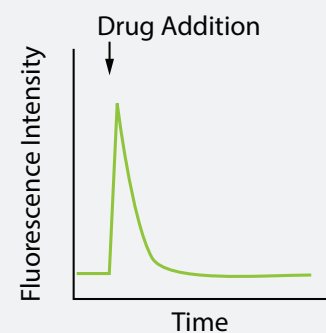
Rq (M1 muscarinic control receptor) 5µL

complete media 43.15µL

total volume per well 50µL

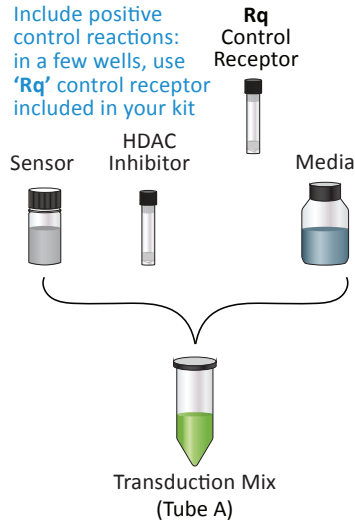
Measure Parameters to Determine Optimal Conditions

- Determine optimal sensor volume by analyzing fluorescence above background, cell health, and response to drug. (For details on measuring fluorescence, refer to [Day 2 procedure](#).)



Day 1 – Transduce and Plate Cells for your Experiment

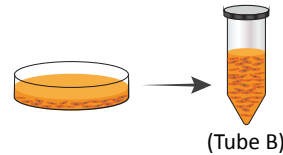
A. **Prepare Viral Transduction Mix (Tube A):** For each transduction reaction (i.e. one well in a 96-well plate), prepare the transduction mix as detailed in table at right (using the optimal volume of sensor that was determined in your optimization experiment). Mix gently.



Transduction Mix (Tube A)	
GE _{CO} Ca ²⁺ Sensor	Variable
SB 500 mM sodium butyrate	0.6μL
Rq control receptor (if using)	5μL
complete media	(To 50μL)
total volume per well	50μL

Tip: When preparing a master transduction mix, to avoid coming up short, scale up by 10-15% of the number of wells needed.

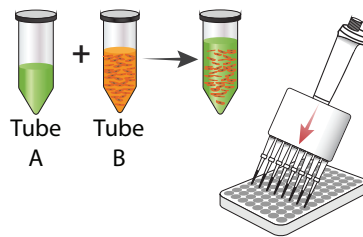
B. **Prepare Cells (Tube B):** Detach cells using standard trypsinization protocol. Resuspend cells in complete culture media and determine cell count.



Cells (Tube B)	
cells per well	100μL

C. **Prepare a dilution of cells at your desired concentration.** (100μL of this cell resuspension will be required for a single well in a 96-well plate, so prepare enough of the dilution to seed the desired number of wells in the plate).

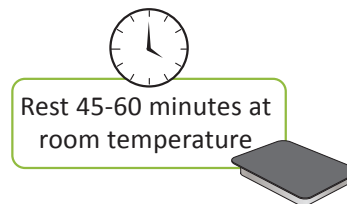
D. **Combine Transduction Mix and Cells:** Combine Tube A and Tube B (50μL Tube A + 100μL Tube B). Mix by pipetting up and down gently, and seed 150μL per well on the 96-well plate.



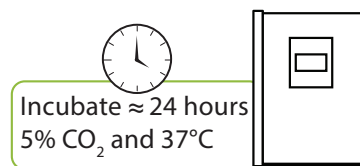
Transduction Mix + Cells (Tube A) + (Tube B)	
transduction mix	50μL
cell suspension	100μL
total volume per well	150μL

Tip: 90% cell confluence after 24 hours is ideal. For HEK293T cells 50,000 cells per well works well but plating density is cell type dependent

E. Cover plate to protect from light and let rest at room temperature for 45-60 minutes.



F. Incubate ≈ 20-24 hours under normal cell growth conditions (5% CO₂ and 37°C), protected from light.





Suggestions for Assays in Adherent Cells

This protocol is optimized for rapidly dividing immortalized cells.

However, GECO Ca²⁺ biosensor assays can be used with primary cultures of cells as well as iPSC-derived cells. In the case of non-dividing, or differentiated cells, the transduction should be done with adherent cells, and the media should be exchanged the following day. Specific details of the protocol will vary by cell type, so it is important to take the time to titrate BacMam for optimal results. For expression in rare cell types, or specific cells in mixed cultures, Cre-dependent and specific promoter systems are available for many of our sensors.

***When working with adherent cell cultures, add an extra day to seed the cells prior to transduction.**

Adherent Cells Day 1:

- Seed the cells the day before you add transduction mix, incubate 24 hours

Adherent Cells Day 2:

- Perform transduction as directed [on page 5](#), but add the transduction reaction directly to the plated cells (no aspiration of cell medium necessary). Gently rock the plate 4-5 times in each direction to mix throughout the well. Incubate the cells under normal growth conditions (5% CO₂ and 37°C), protected from light, for 20-24 hours.
- Optional step (cell type dependent): After 4-8 hour incubation with sensor BacMam (6 hours is optimal), very gently aspirate transduction solution (we recommend using a plate washer). Add 100µL complete growth media with sodium butyrate at a concentration of 2mM. If cells will not tolerate a full media exchange, partial media exchanges can be done.

Adherent Cells Day 3:

- Measure Fluorescence as as detailed in [Day 2 procedure](#).

Scaling for 384 Wells (1 plate)

To set up the assay in a 384-well plate, follow all of the protocol steps, adjusting reagent volumes as follows:

Transduction Mix (Tube A) (384-well plate)

GECO Ca ²⁺ Senso	Variable
500 mM SB sodium butyrate	0.1µL
Rq control receptor	1µL
complete media	(To 12.5µL)

total volume per well 12.5µL

Cells (Tube B) (384-well plate)

cells per well	12.5µL
----------------	--------

12.5 µL of the cell resuspension will be required for a single well in a 384-well plate. A plating density of 7,500 cells per well is a good starting point, so prepare the cell suspension at 600,000 cells/mL. Depending on the cell type and plate type, 5,000-15,000 cells per well may be optimal.

Transduction Mix + Cells (Tube A) + (Tube B) (384-well plate)

transduction mix	12.5µL
cell suspension	12.5µL

total volume per well 25µL

Tip: When scaling for 384-well plates, the volume of the cell suspension per well can be increased to 50µL to improve cell health. Make sure that the well is still receiving 7,500 cells and that you increase the volume of sodium butyrate to 0.2µL per well.



Alternative Spinoculation Protocol

This alternative protocol may also be useful for particularly difficult-to-transduce or low expressing cell types:

- a. Prepare transduction mix (detailed in table at right).
- b. Detach cells using standard trypsinization protocol. Resuspend cells in complete culture media and perform cell count.
- c. Prepare a dilution of cells at your desired concentration (22,500 cells/well in a 96-well plate is a good starting point, but will ultimately depend on the cell type being used). 50µL of this cell resuspension will be required for a single well in a 96-well plate, so prepare enough of the dilution to seed the desired number of wells in the plate.
- d. Combine the transduction mix with the cell suspension (50µL transduction mix + 50µL cells). Mix gently, then seed 100µL of this mix per well on a 96-well plate.
- e. Let cells sit at room temperature, protected from light, for 20 minutes.
- f. **Spin the plate at 1,500 x g for 1.5-2 hours at room temperature.**
- g. * We recommend sealing the plate with Breathe-Easy® (Cat. No. 70536-10) during this step to avoid contamination.
- h. After spinning the plate, **remove the transduction mix** and replace with fresh media containing **0.6µL sodium butyrate** (2mM per well).
- i. Return plate to normal growth conditions and incubate for 48 hours.

Alternative Transduction Mix	
GECO Ca ²⁺ Sensor	Variable
500 mM SB sodium butyrate	0.6µL
1M HEPES (pH 7.4)	0.7µL
cell culture media	(To 50µL)
total volume per well	50µL

Tip: Titrate the sensor to determine optimal volume; see [optimization page](#)

If the above protocol does not result in acceptable expression levels, we recommend using **System Biosciences' Spinoculation protocol**, which adds SBI's TransDux Max and MAX enhancer reagents to the transduction reaction. Please make the following adjustments to your transduction reaction:

Transduction Mix + Enhancer Reagents	
GECO Ca ²⁺ Senso	Variable
TransDux	0.34µL
Max Enhancer	17µL
500 mM SB sodium butyrate)	0.6µL
1M HEPES (pH 7.4)	0.7µL
cell culture media	(To 50µL)
total volume per well	50µL

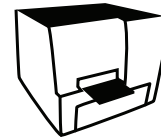
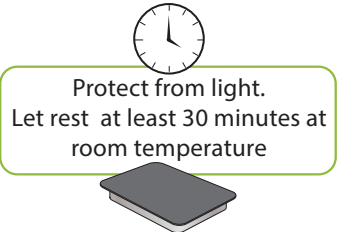
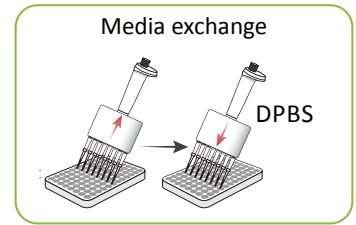


Day 2– Measure Ca²⁺ Response

- G. Prior to measuring fluorescence, replace culture media with DPBS (1X, containing Ca²⁺ and Mg²⁺). Wash gently 4-5 times, careful not to dislodge cells. We recommend using a plate washer. (We recommend the [BioTek 405 TS](#).)
- H. Cover the cells. Allow to equilibrate at room temperature at least 30 minutes before measuring fluorescence. Experiments can be performed at room temperature.
- I. Visually inspect cells on microscope to confirm cell health. Transduced cells should be at least 5× brighter compared to untransduced cells. (If you have high background fluorescence, gently wash cell culture media again.)

Measure Fluorescence on Plate Reader or Imaging System

- J. Measure fluorescence in transduced and untransduced cells. Transduced cells should be at least 5× brighter compared to untransduced cells.
- K. Acquire 10-20 baseline fluorescence reads before adding compounds.
- L. Activate the receptor with agonist to affect levels of Ca²⁺. Resume measurement immediately after adding drug (within 1-2 seconds). Measure continuously for a minimum of 3 minutes. A change in fluorescence intensity under standard GFP excitation and emission will indicate a change in Ca²⁺ levels. For wells transduced with the M1 muscarinic control receptor, add 10μM **Aq** (carbachol, final concentration in well) to activate the receptor. The optimal dose of agonist may need to be determined for a given cell line.



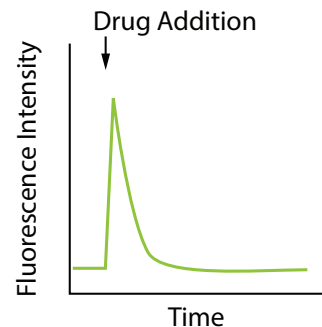
Aq
Control
Agonist



* **Important:** Make sure that the final concentration of any drug solvents in the well (e.g. DMSO) does not exceed 1%. The ideal concentration for DMSO is 0.1%.

Tip: Add agonists at a concentration of 3-4× (in a volume of 50μL DPBS) to wells containing 100-150μL DPBS. This will result in a 1:3 or 1:4 dilution, and allow for adequate diffusion.

Tip: Always test the addition of the vehicle alone (i.e. DPBS without drug).





Fluorescence Detection

Our assays are compatible with automated fluorescent plate readers and imaging systems.

We have validated on:

- Agilent (BioTek) Synergy
- Agilent (BioTek) Cytation™
- BMG CLARIOstar®
- Agilent (BioTek) Lionheart
- Agilent (BioTek) Neo
- Hamamatsu FDSS®
- Epifluorescence microscopes

Our customers have reported good results on:

- Molecular Devices FLIPR®
- Molecular Devices Flexstation®
- Perkin Elmer Enspire®
- Perkin Elmer Opera Phenix®

Tip: Below are the specific settings recommended for use with our assays for instruments we use at Montana Molecular. To determine the best settings for fluorescence detection on your instrument, please consult the manufacturer.

Table 3. CLARIOstar® Recommendations

Instrument Settings	
Detection Mode:	FI (Bottom)
Detection Method:	Plate Mode, Kinetic
Scan Mode:	Orbital Averaging
Scan Diameter (mm):	4
Gain/Focal Height:	Adjusted prior to test
Optical Settings	
Excitation:	F 482-16
Dichroic:	LP 504
Emission:	F 530-40

Table 4. Neo & MX™ Recommendations

Instrument Settings	
Detection Method:	Fluorescence Intensity
Read Type:	Endpoint/Kinetic
Optics:	Monochromators
Excitation:	485 nm
Emission:	528 nm
Bandwidth:	20 nm (for both ex and em)
Optics Position:	Bottom
Gain:	100

Table 5. Cytation™ & Lionheart Recommendations

Image Preprocessing	
Image Set:	GFP
Background:	Dark
Rolling Ball Diameter:	Automatic
Image Smoothing:	0 cycles
Cellular Analysis Parameters	
Channel:	Tsf [GFP]
Threshold:	7,000
Background:	Dark
Split Touching Objects:	Checked
Fill Holes in Mask:	Checked
Minimum Object Size:	5 µm
Maximum Object Size:	1,000 µm
Include Primary Edge Objects:	Checked
Analyze Entire Image:	Checked
Advanced Detection Options	
Rolling Ball Diameter:	Automatic
Image Smoothing Strength:	1 cycle of 3×3 average filter
Evaluate Background On:	5%
Primary Mask:	Use threshold mask



Fluorescence Properties

Green GECO is constructed with a green fluorescent protein. The optimal excitation wavelength for the green fluorescent GECO sensor is 480 nm, but the absorption band of this protein is quite broad, so broad bandpass filters that pass 450 to 480 nm light can be used effectively. On the emission side, the green light spans 510 to 550 nm, so broad band pass emission filters can also collect much of the emission. These filter properties are similar to many of the FITC filter sets commonly available on most microscopes and plate readers. Preferred excitation and emission wavelengths are 488/525..

Tip: We offer [mNeon BacMam Kits](#)

[[products #P1000G \(CMV\)](#), [#P1005G \(CAG\)](#) and [#P1001G \(Synapsin\)](#)]. These test kits are a good way to determine BacMam transduction efficiency, evaluate promoter systems, and optimize expression in your cells of choice.

Timing

Unlike many assays that measure accumulation of Ca^{2+} in cell lysates, the Green GECO assay measures calcium in living cells, in real time. Changes in calcium can occur quite rapidly. For best results, be sure to capture changes in Ca^{2+} during the peak response. In **Figure 2**, fluorescence was captured from cells before the addition of the drug and then sampled at regular intervals. The maximal response is reached at ~2-5 seconds after the addition of the agonist and in HEK 293T cells, returns to baseline after ~60 seconds.

Data Analysis

Check out our helpful reference document "[Biosensor Analysis Overview](#)"

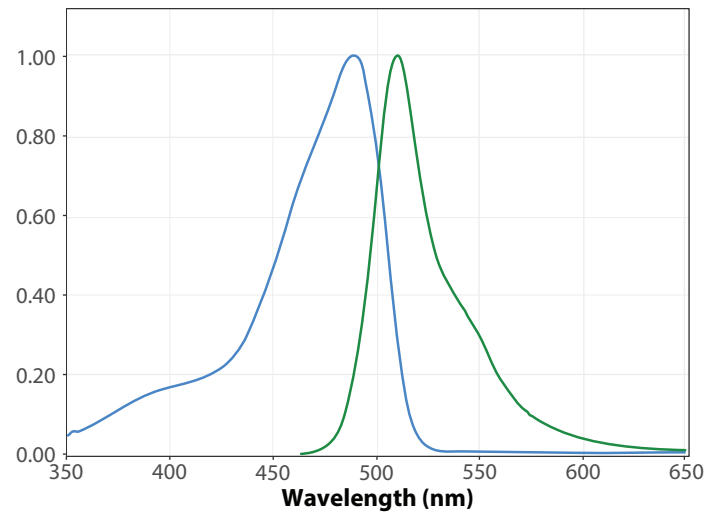


Figure 1. Absorption and emission properties of a typical green fluorescent protein plotted as a function of wavelength.

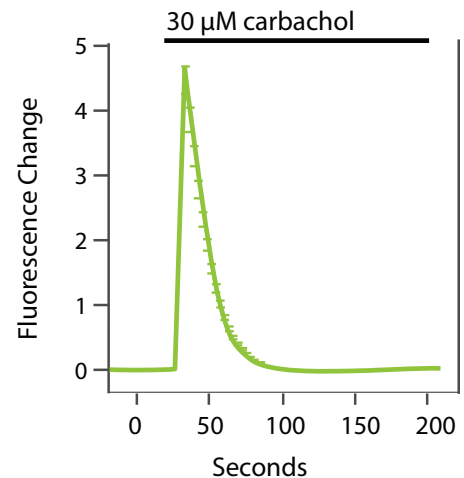


Figure 2. HEK 293T cells transduced with the Green GECO biosensor, and activated with carbachol. The graph shows the expected timing and duration of the Ca^{2+} sensor when using the positive controls provided in your kit.



Assay Optimization

Optimizing Fluorescence

Twenty-four hours after transduction, check your cells for fluorescence. Wells that were transduced with the sensor should be 5-10 times brighter than control wells that were not treated with the sensor.

HDAC inhibitors may be used to maintain expression of the sensors. While BacMam transduction alone will result in sensor expression, sodium butyrate or another HDAC inhibitor, such as valproic acid (VPA) or trichostatin A (TSA), will generate higher levels of expression and will maintain this level of expression [6]. If cells look unhealthy, use lower concentrations or no HDAC inhibitor.

The type of cell culture media used in your experiment can affect the transduction efficiency of BacMam. Our assays have been validated in EMEM, McCoy's 5A, and F12K culture media. Cell culture media with high sodium bicarbonate content can interfere with transduction efficiency [8]. If your cell culture media of choice is affecting transduction efficiency, transduction can be conducted in DPBS for 6 hours and then the cells can be returned to complete cell culture media containing 2mM sodium butyrate.

Optimizing Expression for Your Cell Type

To determine optimal conditions for your cell type, prepare a dilution series of transduction reactions by varying the amount of sensor. For example, a range of 1.25 μ L to 20 μ L is a good starting range (for Product #U0650G) in a 96-well plate. Choose the range that gives you at least 5-10 times above background and gives you the highest amount of fluorescence change with your Rq control receptor.

Varying the cell density, concentration of sodium butyrate, or trying a new HDAC inhibitor (VPA or TSA) may boost expression as well.

Please [contact us](#) if you would like to use the sensor under the control of a specific promoter system. Sensors under weak promoters may be limited to detection on imaging systems. To maintain strong expression in specific cell types, we recommend ordering a Cre-inducible, floxed sensor.

Purified viral preparations, which can increase expression in particularly sensitive or difficult to transduce cell types, are available upon request.

Use the Positive Control

If the cells are expressing the sensor, and fluorescence is detectable on your instrument, then evaluate sensor performance using the positive control receptor included in your kit. Add 5 μ L of the Rq (M1 control receptor) to a set of control wells and activate with isoproterenol included in the kit.

Optimizing Receptor Expression

If you have titrated the GECO sensor and determined the optimal volume, but fail to see a receptor mediated signal, the receptor expression level may be the issue. Try titrating the receptor with a fixed amount of the GECO sensor.

If you need further help, see the [Troubleshooting Guide](#) or let us know, we're happy to help!

+1 406-200-8321, info@montanamolecular.com



Troubleshooting Guide

Problem	Possible Cause	Solution
1. Low sensor expression and/or poor transduction efficiency	Suboptimal sensor BacMam volume is being used.	Perform titration of the sensor BacMam stock, testing a large range (i.e. 2.5-40µL in 96-well plate format) to identify optimal volume. Too little can result in low expression, too much can cause cells to become sick.
	Transducing adherent cells.	Transduce cells while in suspension. If this isn't possible, try doing a media exchange on adherent cells after 4-6 hours, in addition to leaving the virus on overnight. See Suggestions for Assays in Adherent Cells .
	Suboptimal cell density; too few or too many cells added.	Transduce cells so that the cells will be around 75-80% confluent at the time of transduction. Also, when transducing cells in suspension, make sure that cells in the source flask are < 100% confluent (approximately 80% confluent is ideal).
	HDAC inhibitor was not added to the transduction mix, or the concentration was wrong.	Add HDAC inhibitor at the proper concentration: sodium butyrate - 2 mM valproic acid - 5 mM trichostatin A - 0.25µM * Perform a titration to determine optimal concentration for the cell type being used.
	HDAC inhibitor being used is not optimal for cell type.	Test other HDAC inhibitors (e.g. sodium butyrate, valproic acid, trichostatin A.)
	Cell type being used transduces poorly.	<ul style="list-style-type: none"> After adding transduction mix to cells, let cells sit at room temperature for 30-40 min. before placing back in incubator (longer incubation times at room temperature may further improve transduction). Perform media exchange after various incubation times with the transduction mix, in addition to leaving the virus on overnight. Try high-titer, purified BacMam stock. Validate assay in a different cell type (e.g. HEK 293T cells) Transduce cells multiple times (e.g. on Day 1, and again on Day 2). Incubate cells for 48 hours post transduction, before performing assay. Consider using a different viral vector, such as lentivirus or AAV.
	Cell culture media is inhibiting transduction.	Remove media during transduction, preparing the transduction mix in PBS and adding to cells. Replace transduction mix with media after 2-4 hours.
	BacMam stock was not stored properly (i.e. not stored at 4°C, exposed to light for long periods, subjected to multiple freeze-thaw cycles), or the shelf life has been exceeded.	Follow guidelines for product storage . BacMam stocks are stable for at least 12 months when stored properly. After this time period, the stock should be re-evaluated and compared to previous experiments. Purified BacMam stocks should be used within 30 days for best results.
	BacMam stock was not mixed adequately before transducing cells	Mix BacMam stock thoroughly before transduction, especially after being stored for long periods.
	Promoter is not optimal for cell type being used	Identify promoters that work best in the cell type being used. If promoter is not on product list, consult Montana Molecular for custom production services.
Cells are contaminated	Monitor cells for bacteria, fungi, etc	



Problem	Possible Cause	Solution
2. Low fluorescence signal on microscope/plate reader.	Low sensor expression, low transduction efficiency.	See solutions for Problem 1 .
	Excitation/emission settings are not optimal for the sensor being used.	Refer to protocol for the fluorescence spectra of the sensor. Make sure that filter sets or monochromators are aligned with the peak excitation and emission wavelengths of the sensor.
	Cells are in cell culture media, and the media is producing a large fluorescent signal (autofluorescence).	Exchange media so that cells are in PBS at the time of experiment. If cell culture media must be used, try using FluoroBrite media.
	Wrong microplate type is being used.	Use black, clear-bottom microplates with low autofluorescence.
	Exposure time or gain setting on instrument is suboptimal.	Test different exposure and gain settings, monitoring how the signal-to-background and noise in the measurement changes. Too high can result in saturation and/or photobleaching; too low can result in noisy data.
	Cells were dislodged during media exchange/plate washing.	Make sure that media exchange or plate washing is done gently and does not dislodge cells. Confirm with visual inspection on a microscope.
3. Signal-to-background is low (i.e. cells/wells with sensor are not much brighter than control cells/wells without sensor).	Low sensor expression, low transduction efficiency.	See solutions for Problem 1 .
	Excitation/emission settings are not optimal for the sensor being used.	Refer to protocol for the fluorescence spectra of the sensor. Make sure that filter sets or monochromators are aligned with the peak excitation and emission wavelengths of the sensor.
	Exposure time or gain setting on instrument is suboptimal.	Test different exposure and gain settings, monitoring how the signal-to-background and noise in the measurement changes. Too high can result in saturation and/or photobleaching; too low can result in noisy data.
	Media exchange was not performed before running the assay; cells are in media rather than PBS. Cell culture media being used has high autofluorescence.	Perform media exchange so that cells are in PBS at the time of experiment. If cell culture media must be used, try using FluoroBrite media.
	Cells were dislodged during media exchange/plate washing.	Make sure that media exchange or plate washing is done gently and does not dislodge cells. Confirm with visual inspection on a microscope.



Problem	Possible Cause	Solution
4. Signal is noisy.	Low sensor expression, low transduction efficiency.	See solutions for Problem 1 .
	Gain setting or exposure time on instrument is too low.	Increase gain setting or exposure time.
	Media exchange was not performed, or plate washing was inadequate causing high well-to-well variability. Cells are not in PBS at the time of experiment.	Exchange media so that cells are in PBS at the time of experiment. If cell culture media must be used, try using FluoroBrite media. Make sure that plate washing is highly consistent from well to well.
	Cells were dislodged during media exchange/plate washing.	Make sure that media exchange or plate washing is done gently and does not dislodge cells. Confirm with visual inspection on a microscope.
5. Good fluorescence signal, but sensor is not responding to drug as expected. No change in fluorescence observed, or signal is in the wrong direction.	Photobleaching	Reduce exposure time, sampling rate, and/or light intensity.
	Drug is at the wrong concentration	Confirm drug concentration and solubility.
	Drug was not stored properly.	Confirm drug storage conditions.
	Drug was added to the cells in a volume that was too low relative to the volume of PBS/media in the well, resulting in improper mixing.	Add drug in a volume that will allow for sufficient diffusion (i.e. 1:3 or 1:4 drug to total volume)
	Drug was not added in the same solution as the solution in the well/culture dish.	Make sure that the drug preparation and cells are in the same solution.
	Drug addition is producing an artifact.	Make sure to add a vehicle-only control. Make sure drug is added in a solution that is the same as the solution in the well. Do not exceed 1% DMSO final in the well (0.5% or less is ideal).
	Compounds being tested are fluorescent.	Scan compounds for fluorescence to confirm. If possible, dilute compounds in order to reduce the fluorescence artifact of the compound.
	Drug addition was too forceful and dislodged cells.	Add drugs manually or with an on-board dispense function, but do so gently, so as not to dislodge cells.
	Baseline reads were not acquired before adding drug.	Acquire 5-10 baseline fluorescence reads before adding drug. Monitor for a change in fluorescence intensity upon addition of drug.
	Gain setting on instrument is too high, and signal is saturating. Gain setting is too low, and signal cannot be detected.	Adjust gain setting.
	Too much sensor has been added to cells and the signal is saturated (i.e. not enough analyte for the amount of sensor in the cell).	Titrate the amount of sensor to determine maximum signal for your cell type. See protocol recommendations for HEK 293T cells .
	Target receptor was not added, or expression levels are suboptimal (too little or too much, or receptor has high level of constitutive activity).	Titrate the amount of receptor to optimize the signal for your cell type and receptor.
Sampling rate is not consistent with sensor kinetics.	Acquire 5-10 baseline measurements before adding drug. Resume measurement quickly after adding drug (within 5-10 seconds for DAG/PIP ₂ , 60 seconds for cADDi and cGMP, and 1-2 seconds for GECO Ca ²⁺). Measure long enough to capture max response of sensor.	



Problem	Possible Cause	Solution
6. Poor cell health, cells detaching from plate.	Too much BacMam stock was added to cells (e.g. sensor, receptor, Gs mutant).	Titrate lower amounts of BacMam stock to identify the optimal volume for your cells.
	Concentration of HDAC inhibitor is too high, or cells are sensitive to the HDAC inhibitor being used.	Confirm concentration of HDAC inhibitor being used. Make new stock solution. Try a different HDAC inhibitor. Confirm that they are being used at the proper concentration: sodium butyrate - 2 mM valproic acid - 5 mM trichostatin A - 0.25 μM * Perform a titration to determine optimal concentration for the cell type being used.
	Plate surface is not coated with a cell attachment factor.	Coat plates with a cell attachment factor (e.g. PDL, laminin, collagen, fibronectin etc.) to enhance attachment.
	Edge wells are being used, and cells in the edge wells may be subject to conditions that are not conducive to growth.	Do not use edge wells.
	Cells were dislodged during media exchange/plate washing.	Make sure that media exchange or plate washing is done gently and does not dislodge cells. Confirm with visual inspection on a microscope.
	PBS being used does not contain Ca ²⁺ and Mg ²⁺ .	Use PBS containing Ca ²⁺ and Mg ²⁺ .
	Cells are contaminated.	Monitor cells for bacteria, fungi, mycoplasma.
	Cells were not grown under proper growth conditions (i.e. 5% CO ₂ , 37°C).	Incubate transduced cells at 37°C, in 5% CO ₂ .
Cells are sensitive to acidity from the insect cell culture media present in BacMam virus. The insect cell culture media is more acidic than typical mammalian cell culture media.	Wash transduction mix off of cells after 4-6 hours. Replace with fresh cell culture media, maintaining concentration of sodium butyrate or other HDAC inhibitor.	



GECO Calcium Sensor - in the Literature

1. M.C. Cañizal, et al. [A Dual Ca²⁺/ DAG Sensor Reports on Ligand Efficiency](#).
2. C. Xingjuan, et al. [PKC-dependent Phosphorylation of the H1 Histamine Receptor Modulates TRPC6 Activity](#). Cells. 2014.
3. P. Tewson, et al. [A multiplexed fluorescent assay for independent second-messenger systems: decoding GPCR activation in living cells](#). Journal of Biomolecular Screening 18, 2013.
4. J. Wu, et al. [Improved Orange and Red Ca²⁺ Indicators and Photophysical Considerations for Optogenetic Applications](#). ACS Chem Neurosci. Jun. 19, 2013
5. H.J. Carlson, et al. [Mutational Analysis of a Red Fluorescent Protein-Based Calcium Ion Indicator](#). Sensors, 2013.
6. P. Tewson, et al. [Simultaneous Detection of Ca²⁺ and Diacylglycerol Signaling in Living Cells](#). PLoS One. 2012.

References

1. Zhao Y, Araki S, Wu J, et al. [An Expanded Palette of Genetically Encoded Ca²⁺ Indicators](#). Science 2011; 333(6051):1888-1891. doi:10.1126
2. Boyce, FM; Bucher, NL (1996). [Baculovirus-mediated gene transfer into mammalian cells](#). Proceedings of the National Academy of Sciences of the United States of America 93 (6): 2348–52. doi:10.1073/pnas. 93.6.2348. PMC 39799. PMID 8637876.
3. Graham FL, Smiley J, Russell WC, Nairn R: [Characteristics of a human cell line transformed by DNA from human adenovirus type 5](#). J Gen Virol 1977, 36(1):59-74.
4. Dulbecco R and Vogt M: [Plaque formation and isolation of pure lines with poliomyelitis viruses](#). The Journal of experimental medicine 1954.
5. Tewson, et al. [Simultaneous Detection of Ca²⁺ and Diacylglycerol Signaling in Living Cells](#). PLoS One, 2012.
6. Kost T, Condreay J, Ames R, Rees S, Romanos M: [Implementation of BacMam virus gene delivery technology in a drug discovery setting](#). Drug Discovery Today 2007, 12(9-10):396-403
7. Pidre, M. L., Arrías, P. N., Amorós Morales, L. C., & Romanowski, V. (2022). The Magic Staff: A Comprehensive Overview of Baculovirus-Based Technologies Applied to Human and Animal Health. Viruses, 15(1). <https://doi.org/10.3390/v15010080>
8. Shen H-C, et al., [Baculovirus-mediated gene transfer is attenuated by sodium bicarbonate](#). J Gene Med. Jun. 2007.

Related Products

Product	Sensor Description	Promoter	Recommended Use
#U0250R	Red Up cADDIs cAMP - Big Sky	CMV	Fluorescence imaging and plate read assay
#D0350R	Red Down DAG - Big Sky	CMV	Fluorescence imaging and plate reader assay
#D0450R	Red Down PIP2 Sensor - Big Sky	CMV	Fluorescence imaging

Contact Us

If you have any questions about the protocols described here, or if you have ideas about how we can improve these tools, then we want to hear from you. Your feedback is extremely valuable. Please send an email to info@montanamolecular.com or call us at +1 406-200-8321 and we'll respond as quickly as we can.

