



Product Code	Fluorescent Biosensor BacMam Kit	Promoter	Use	Sizes	List Price (USD)
#D0200G #D0205G	Green Downward cADDis cAMP	CMV CAG	Screen Gs-coupled GPCRs to monitor increases in cAMP on fluorescence plate readers or imaging systems	10mL 30 mL 60 mL	\$595 \$1195 \$1795
#U0200G #U0205G	Green Upward cADDis cAMP	CMV CAG	Screen Gs-coupled GPCRs to monitor increases in cAMP on fluorescence plate readers or imaging systems	10mL 30 mL 60 mL	\$595 \$1195 \$1795
#U0200R #U0205R	Red Upward cADDis cAMP	CMV CAG	Screen Gs-coupled GPCRs to monitor increases in cAMP on fluorescence plate readers or imaging systems	10mL 30 mL 60 mL	\$595 \$1195 \$1795
#X0200G #X0205G	Green Gi cADDis cAMP	CMV CAG	Screen Gi-coupled GPCRs to monitor cAMP reduction on fluorescence plate readers or imaging systems	10mL 30 mL 60 mL	\$595 \$1195 \$1795
#D0201G	Green Cilia-Targeted cADDis cAMP	CMV	Monitor changes in cAMP in the primary cilia with high N.A. fluorescence microscopy.	10mL 30 mL	\$595 \$1195
#D0211G	Ratiometric Cilia-Targeted cADDis cAMP	CMV	Monitor changes in cAMP in the primary cilia with high N.A. fluorescence microscopy.	10mL 30 mL	\$595 \$1195
#D0300G #D0305G	Green Downward DAG	CMV CAG	Screen Gq-coupled GPCRs to monitor increases in DAG on fluorescence plate readers or imaging systems	10mL 30 mL 60 mL	\$595 \$1195 \$1795
#U0300G #U0305G	Green Upward DAG	CMV CAG	Screen Gq-coupled GPCRs to monitor increases in DAG on fluorescence plate readers or imaging systems	10mL 30 mL 60 mL	\$595 \$1195 \$1795
#D0308G	Cre-Dependent Green Downward DAG	Ef1- α	Target expression to Cre-expressing cells to monitor diacylglycerol on fluorescence imaging systems	10mL 30 mL	\$595 \$1195
#D0300R	Red Downward DAG	CMV	Monitor Gq-mediated changes in diacylglycerol on fluorescence imaging systems	10mL 30 mL 60 mL	\$595 \$1195 \$1795
#U0300R	Red Upward DAG	CMV	Monitor Gq-mediated changes in diacylglycerol on fluorescence imaging systems	10mL 30 mL	\$595 \$1195
#D0400G #D0405G	Green PIP ₂ Sensor	CMV CAG	Monitor changes in PIP ₂ on fluorescence imaging systems	10 mL	\$595
#D0400R	Red PIP ₂ Sensor	CMV	Monitor changes in PIP ₂ on fluorescence imaging systems	10 mL	\$595
#C0500N #C0505N #C0506N	Cre Recombinase	CMV CAG Synapsin	Expresses Cre	10 mL	\$395

Specialized promoter systems, lenti and AAV vectors are [available by request.](#)

Product Code	Fluorescent Biosensor BacMam Kit	Promoter	Use	Sizes	List Price (USD)
#U0600R	Red GECO Ca ²⁺	CMV	Red fluorescent Ca ²⁺ indicator can be paired with green cADDIs or DAG to detect simultaneous second messenger signals.	10mL 30 mL	\$595 \$1195
#U0608G	Green GECO Ca ²⁺	Ef1- α	Green Ca ²⁺ indicator with Ef1- α promoter.	10mL 30 mL	\$595 \$1195
#D0700G	ArcLight Voltage	CMV	Monitor electrical activity in excitable cells with changes in green fluorescence.	10 mL	\$595
#D0800G	Green Downward GENle cGMP	CMV	Detect changes in cGMP with a genetically encoded nucleotide indicator (GENle).	10 mL	\$595
#U0900G #U0901G	Green Cell Stress Green Cell Stress Red/Green Ratiometric	CMV	Bright fluorescent assay to detect chemical or genetic ER stress response and UPR.	10mL 30mL 60mL	\$595 \$1195 \$1795

Product Code	GPCR Receptors in BacMam	Promoter	G-Protein Coupled	Sizes	List Price (USD)
#Z1000N	D2 Dopamine Receptor	CMV	Gi	5mL	\$295
#Z1005N	D2 Dopamine Receptor	CAG	Gi	5mL	\$295
#Z2000N	Human M2 Muscarinic Receptor	CMV	Gi	5mL	\$295
#Z2005N	Human M2 Muscarinic Receptor	CAG	Gi	5mL	\$295
#Z3000N	A1 Adenosine Receptor	CMV	Gi	5mL	\$295
#Z4000N	Succinate Receptor	CMV	Gi	5mL	\$295
#Z5000N	β 2 Adrenergic Receptor	CMV	Gs	5mL	\$295
#Z5005N	β 2 Adrenergic Receptor	CAG	Gs	5mL	\$295
#Z6000N	Human M1 Muscarinic Receptor	CMV	Gq	5mL	\$295
#Z7000N	D1 Dopamine Receptor	CMV	Gs/a	5mL	\$295
#Z8000N	Calcitonin Receptor	CMV	Gs/q	5mL	\$295

Specialized promoter systems, lenti and AAV vectors are [available by request](#).