

Product Code	Promoter	Arrestin Recruitment Biosensors	Use	Sizes	List Price (USD)
#D2000G	CMV	Borealis – V2R	Kinetic arrestin signaling assay through activation of Vasopressin. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2010G	CMV	Borealis – AT1R	Kinetic arrestin signaling assay through activation of Angiotensin. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2020G	CMV	Borealis – GLP-1R	Kinetic arrestin signaling assay through activation of GLP-1R. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2030G	CMV	Borealis – MOR	Kinetic arrestin signaling assay through activation of μ -Opioid receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2040G	CMV	Borealis – β 2AR	Kinetic arrestin signaling assay through activation of β 2-Adrenergic receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2050G	CMV	Borealis – D1R	Kinetic β -arrestin signaling assay through activation of D ₁ Dopamine receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2060G	CMV	Borealis – OTR	Kinetic arrestin signaling assay through activation of Oxytocin receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2070G	CMV	Borealis – 5-HT2AR	Kinetic arrestin signaling assay through activation of Serotonin 5-HT2A receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2080G	CMV	Borealis – KOR	Kinetic arrestin signaling assay through activation of κ -Opioid receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2090G	CMV	Borealis – PAR2	Kinetic arrestin signaling assay through activation of the PAR2 receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2100G	CMV	Borealis – DOR	Kinetic arrestin signaling assay through activation of the DOR receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2110G	CMV	Borealis – LPA1R	Kinetic arrestin signaling assay through activation of the LPA1 receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2120G	CMV	Borealis – PTH1R	Kinetic arrestin signaling assay through activation of the PTH1 receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2130G	CMV	Borealis – 5-HT2CR	Kinetic arrestin signaling assay through activation of the 5-HT2C receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2140G	CMV	Borealis – GCGR	Kinetic arrestin signaling assay through activation of the Glucagon receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2150G	CMV	Borealis – MC4R	Kinetic arrestin signaling assay through activation of the Melanocortin 4 receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2160G	CMV	Borealis – CCR5	Kinetic arrestin signaling assay through activation of the C-C Chemokine Receptor Type 5. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2170G	CMV	Borealis – PAR1	Kinetic arrestin signaling assay through activation of the Protease-activated Receptor 1. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D2180G	CMV	Borealis – Apelin (APJ/APLNR)	Kinetic arrestin signaling assay through activation of the Apelin Receptor (APJ/APLNR). Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895

Product Code	Promoter	Cell Signaling Biosensors	Use	Sizes	List Price (USD)
#D0200G	CMV	Green Down cADDIs cAMP	Screen Gs-coupled GPCRs to monitor increases in cAMP on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
				200 mL	A.Q.
#D0250G #D0251G	CMV Synapsin	Green Down cADDIs cAMP - Big Sky BacMam	Screen Gs-coupled GPCRs to monitor increases in cAMP on fluorescence plate readers or imaging systems. Packaged in our new vector yielding significantly improved transduction efficiency.	5 mL	\$795
				15 mL	\$1695
				30 mL	\$2895
#U0200G	CMV	Green Up cADDIs cAMP	Screen Gs-coupled GPCRs to monitor increases in cAMP on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
				200 mL	A.Q.
#U0250G #U0251G	CMV Synapsin	Green Up cADDIs cAMP - Big Sky BacMam	Screen Gs-coupled GPCRs to monitor increases in cAMP on fluorescence plate readers or imaging systems. Packaged in our new vector yielding significantly improved transduction efficiency.	5 mL	\$795
				15 mL	\$1695
				30 mL	\$2895
#U0200R	CMV	Red Up cADDIs cAMP	Screen Gs-coupled GPCRs to monitor increases in cAMP on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#U0250R	CMV	Red Up cADDIs cAMP - Big Sky BacMam	Screen Gs-coupled GPCRs to monitor increases in cAMP on fluorescence plate readers or imaging systems. Packaged in our new vector yielding significantly improved transduction efficiency.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#X0200G	CMV	Green Gi cADDIs cAMP	Screen Gi-coupled GPCRs to monitor cAMP reduction on fluorescence plate readers or imaging systems. Packaged in our new vector yielding significantly improved transduction efficiency.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
				200 mL	A.Q.
#X0250G	CMV	Green Gi cADDIs cAMP - Big Sky BacMam	Screen Gi-coupled GPCRs to monitor cAMP reduction on fluorescence plate readers or imaging systems. Packaged in our new vector yielding significantly improved transduction efficiency.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D0201G	CMV	Green Cilia-Targeted cADDIs cAMP	Monitor changes in cAMP in the primary cilia with high N.A. fluorescence microscopy.	10 mL	\$795
				30 mL	\$1695
#D0211G	CMV	Ratiometric Cilia-Targeted cADDIs cAMP	Monitor changes in cAMP in the primary cilia with high N.A. fluorescence microscopy.	10 mL	\$795
				30 mL	\$1695
#D0221G	CMV	Green Nucleus-Targeted cADDIs cAMP	Monitor changes in cAMP in the cell nucleus with fluorescence imaging.	10 mL	\$795
				30 mL	\$1695
#U0221R	CMV	Red Nucleus-Targeted cADDIs cAMP	Monitor changes in cAMP in the cell nucleus with fluorescence imaging.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D0231G	CMV	Green Membrane-Targeted fMP cADDIs cAMP	Flag tagged MyrPalm cADDIs targeted to the plasma membrane.	10 mL	\$795
				30 mL	\$1695
#U0241R	CMV	Red Membrane-Targeted fS15 cADDIs cAMP	Flag tagged S15 cADDIs targeted to the plasma membrane (red upward).	10 mL	\$795
				30 mL	\$1695
#D0241G	CMV	Green Membrane-Targeted fS15 cADDIs cAMP	Flag tagged S15 cADDIs targeted to the plasma membrane (green downward).	10 mL	\$795
				30 mL	\$1695
#U0261G	CMV	AKAP-targeted cADDIs Sensors - AKAP18a	Measure changes in signalsomes, with cADDIs sensor attached directly to different AKAP proteins.	5 mL	\$795
#U0262G	CMV	AKAP-targeted cADDIs Sensors - AKAP18b	Measure changes in signalsomes, with cADDIs sensor attached directly to different AKAP proteins.	5 mL	\$795
#U0263G	CMV	AKAP-targeted cADDIs Sensors - AKAP1	Measure changes in signalsomes, with cADDIs sensor attached directly to different AKAP proteins.	5 mL	\$795
#U0264G	CMV	AKAP-targeted cADDIs Sensors - AKAP79	Measure changes in signalsomes, with cADDIs sensor attached directly to different AKAP proteins.	5 mL	\$795
#U0266G	CMV	AKAP-targeted cADDIs Sensors - AKAP12	Measure changes in signalsomes, with cADDIs sensor attached directly to different AKAP proteins.	5 mL	\$795
#U0267G	CMV	AKAP-targeted cADDIs Sensors - smAKAP	Measure changes in signalsomes, with cADDIs sensor attached directly to different AKAP proteins.	5 mL	\$795



Product Code	Promoter	Cell Signaling Biosensors	Use	Sizes	List Price (USD)
#D0300G	CMV	Green Down DAG	Screen Gq-coupled GPCRs to monitor increases in DAG on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
				200 mL	A.Q.
#D0350G	CMV	Green Down DAG - Big Sky BacMam	Screen Gq-coupled GPCRs to monitor increases in DAG on fluorescence plate readers or imaging systems. Packaged in our new vector yielding significantly improved transduction efficiency.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#U0300G	CMV	Green Up DAG	Screen Gq-coupled GPCRs to monitor increases in DAG on fluorescence plate readers or imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#U0350G	CMV	Green Up DAG - Big Sky BacMam	Screen Gq-coupled GPCRs to monitor increases in DAG on fluorescence plate readers or imaging systems. Packaged in our new vector yielding significantly improved transduction efficiency.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D0308G	Ef1- α	Cre-Dependent Green Down DAG	Target expression to Cre-expressing cells to monitor diacylglycerol on fluorescence imaging systems.	10 mL	\$795
				30 mL	\$1695
#D0300R	CMV	Red Down DAG	Monitor Gq-mediated changes in diacylglycerol on fluorescence imaging systems.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#U0300R	CMV	Red Up DAG	Monitor Gq-mediated changes in diacylglycerol on fluorescence imaging systems	10 mL	\$795
				30 mL	\$1695
#D0331G	CMV	Membrane-Targeted DAG	DAG sensor targeted (CAAX) to the plasma.	10 mL	\$795
				30 mL	\$1695
#D0400G	CMV	Green PIP ₂ Sensor	Monitor changes in PIP ₂ on fluorescence imaging systems	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#D0400R	CMV	Red PIP ₂ Sensor	Monitor changes in PIP ₂ on fluorescence imaging systems	10 mL	\$795
#U0600G #U0608G	CMV Ef1- α	Green GECO Ca ²⁺	Green Ca ²⁺ indicator with CMV or Ef1- α promoter.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
				200 mL	A.Q.
#U0650G	CMV	Green GECO Ca ²⁺ - Big Sky BacMam	Green Ca ²⁺ indicator packaged in our new vector yielding significantly improved transduction efficiency.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#U0600R	CMV	Red GECO Ca ²⁺	Red fluorescent Ca ²⁺ indicator can be paired with green cADDis or DAG to detect simultaneous second messenger signals.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
				200 mL	A.Q.
#U0650R	CMV	Red GECO Ca ²⁺ - Big Sky BacMam	Red fluorescent Ca ²⁺ indicator can be paired with green cADDis or DAG to detect simultaneous second messenger signals. Packaged in our new vector yielding significantly improved transduction efficiency.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
#U0651R	CMV	Red Mitochondria-Targeted GECO Ca ²⁺	Red fluorescent Ca ²⁺ indicator, targeted to the mitochondria, can be paired with green cADDis or DAG to detect simultaneous second messenger signals. Packaged in our new vector yielding significantly improved transduction efficiency.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
				200 mL	A.Q.
#D0800G	CMV	Green Down GENIE cGMP	Detect changes in cGMP with a genetically encoded nucleotide indicator (GENIE).	30 mL	\$1695

Product Code	Promoter	Cell Stress Biosensors	Use	Sizes	List Price (USD)
#U0900G	CMV	Green Cell Stress	Bright fluorescent assay to detect response and reversal of chemical or genetic ER stress and UPR.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
				200 mL	A.Q.
#U0901G	CMV	Green/Red Ratiometric Cell Stress	Bright fluorescent assay to detect response and reversal of chemical or genetic ER stress and UPR.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
				200 mL	A.Q.
#U0921G	CMV	IRE1-XBP1 Cell Stress	Bright fluorescent assay to detect response and reversal of chemical or genetic ER stress and UPR.	10 mL	\$795
				30 mL	\$1695
				60 mL	\$2895
				200 mL	A.Q.

Product Code	Promoter	SARS-CoV-2 Reagents	Use	Sizes	List Price (USD)
#C1160G	CMV	3CLglow	Green fluorescent biosensor for the SARS-CoV-2 Main protease. Screen for potential inhibitors of SARS-CoV-2 in living cells.	5 mL 30 mL 60 mL	\$695 \$2595 \$3895
#C1161G	CMV	3CLglowUp	Our newest and brightest tool for screening 3CL protease inhibitors. 3CLglowUp sensor increases in fluorescence in response to inhibition of the 3CL ^{pro} enzyme.	5 mL 30 mL 60 mL	\$695 \$2595 \$3895
#C1162G	CMV	3CLglowLive	The 3CLglowLive Assay kit is well suited to experiments using the live SARS-CoV-2 virus. It is dim at baseline, but fluoresces green in the presence of 3CL protease activity, an indicator of SARS-CoV-2 viral replication.	5 mL 30 mL 60 mL	\$695 \$2595 \$3895
#C1100G	CMV	mNeonGreen Fluorescent ACE2	Controlled expression of green fluorescent ACE2 for SARS-CoV-2 drug screening.	5 mL 30 mL 60 mL 200 mL	\$395 \$795 \$1195 A.Q.
#C1100R	CMV	Red Fluorescent ACE2	Controlled expression of red fluorescent ACE2 for SARS-CoV-2 drug screening.	5 mL 30 mL 60 mL 200 mL	\$395 \$795 \$1195 A.Q.
#C1100N	CMV	Unlabeled ACE2	Controlled expression of (unlabeled) ACE2 for SARS-CoV-2 drug screening.	5 mL 30 mL 60 mL 200 mL	\$395 \$795 \$1195 A.Q.
#C1110G	CMV	Pseudo SARS-CoV-2 Green Reporter	BacMam virus pseudotyped with SARS-CoV-2 spike protein that will produce green fluorescent nuclei upon successful entry into host cells expressing ACE2.	5 mL 30 mL 60 mL 200 mL	\$695 \$2595 \$3895 A.Q.
#C1120G	CMV	Pseudo SARS-CoV-2 D614G Green Reporter	BacMam virus pseudotyped with SARS-CoV-2 D614G mutated spike protein that binds with ACE2 and expresses green fluorescent protein at the nucleus 24 hours after viral entry.	5 mL 30 mL 60 mL 200 mL	\$695 \$2595 \$3895 A.Q.
#C1121G	CMV	Pseudo SARS-CoV-2 Spike M1 Green Reporter	BacMam virus pseudotyped with SARS-CoV-2 D614G, E484K, N501Y, K417T mutated spike protein that binds with ACE2 and expresses green fluorescent protein at the nucleus 24 hours after viral entry.	5 mL 30 mL 60 mL 200 mL	\$695 \$2595 \$3895 A.Q.
#C1122G	CMV	Pseudo SARS-CoV-2 Spike M2 Green Reporter	BacMam virus pseudotyped with SARS-CoV-2 D614G, E484K, N501Y, K417N mutated spike protein that binds with ACE2 and expresses green fluorescent protein at the nucleus 24 hours after viral entry.	5 mL 30 mL 60 mL 200 mL	\$695 \$2595 \$3895 A.Q.
#C1123G	CMV	Pseudo SARS-CoV-2 Green Reporter with Spike Delta Variant	The Delta Variant of the Spike protein, is also known as B.1.617.2 and includes the following Spike protein mutations: T19R, V70F, T95I, G142D, E156-, F157-, R158G, A222V, W258L, K417N, L452R, T478K, D614G, P681R, D950N.	5 mL 30 mL 60 mL 200 mL	\$695 \$2595 \$3895 A.Q.
#C1130N	CMV	TMPRSS2	TMPRSS2 in BacMam for controlled expression in live mammalian cells and increased spike binding.	5 mL 30 mL 60 mL	\$395 \$795 \$1195
#C1140R	CMV	Neuropilin-1 - Red	Red fluorescent labeled Neuropilin 1 in a BacMam vector for controlled expression in a wide variety of cell types	5 mL 30 mL	\$395 \$795
#C1140N	CMV	Neuropilin-1 - Unlabeled	Unlabeled Neuropilin 1 in a BacMam vector for controlled expression in a wide variety of cell types.	5 mL 30 mL	\$395 \$795
#C1150N	CMV	Cathepsin-L	Cathepsin L in BacMam for efficient delivery to practically any cell type.	5 mL 30 mL	\$395 \$795
#C1200N	CMV	3CL ^{pro} T21I Expression Kit	Express 3CL ^{pro} with the T21I mutation in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1201N	CMV	3CL ^{pro} L50F Expression Kit	Express 3CL ^{pro} with the L50F mutation in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1202N	CMV	3CL ^{pro} E166V Expression Kit	Express 3CL ^{pro} with the E166V mutation in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1203N	CMV	3CL ^{pro} E166G Expression Kit	Express 3CL ^{pro} with the E166G mutation in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1204N	CMV	3CL ^{pro} Q189E Expression Kit	Express 3CL ^{pro} with the Q189E mutation in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1205N	CMV	3CL ^{pro} H172Y Expression Kit	Express 3CL ^{pro} with the H172Y mutation in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1206N	CMV	3CL ^{pro} E166A + L167F Expression Kit	Express 3CL ^{pro} with the E166A & L167F mutations in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1207N	CMV	3CL ^{pro} T21I + T304I Expression Kit	Express 3CL ^{pro} with the T21I & T304I mutations in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1208N	CMV	3CL ^{pro} T21I + E166V Expression Kit	Express 3CL ^{pro} with the T21I & E166V mutations in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1209N	CMV	3CL ^{pro} L50F + E166V Expression Kit	Express 3CL ^{pro} with the L50F & E166V mutations in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1210N	CMV	3CL ^{pro} L50F + E166A + L167F Expression Kit	Express 3CL ^{pro} with the L50F, E166A & L167F mutations in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1211N	CMV	3CL ^{pro} T21I + S144A + T304I Expression Kit	Express 3CL ^{pro} with the T21I, S144A & T304I mutations in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1212N	CMV	3CL ^{pro} T21I + A173V + T304I Expression Kit	Express 3CL ^{pro} with the T21I, A173V & T304I mutations in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1213N	CMV	3CL ^{pro} T21I + L50F + A193P + S301P Expression Kit	Express 3CL ^{pro} with the T21I, L50F, A193P & S301P mutations in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1214N	CMV	3CL ^{pro} L50F + F140L + L167F + T304I Expression Kit	Express 3CL ^{pro} with the L50F, F140L, L167F & T304I mutations in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795
#C1215N	CMV	3CL ^{pro} T21I + C160F + A173V + V186A + T304I Expression Kit	Express 3CL ^{pro} with the T21I, C160F, A173V, V186A & T304I mutations in mammalian cells. Use with other mutant proteases or 3CLglow biosensor to test 3CL ^{pro} inhibitor efficacy against a panel of mutations in the protease.	5 mL 30 mL	\$395 \$795



Product Code	Promoter	Cell Painting Fluorescent Proteins	Use	Sizes	List Price (USD)
#P1000R	CMV	Untargeted Montana Paintbrush (MTPaint)	A quickly folding bright cytosolic red fluorescent protein in a BacMam vector for controlled expression in a wide variety of cell types.	5 mL	\$595
#P1010R	CMV	Mitochondrial-targeted Montana Paintbrush (MitoPaint)	The inner mitochondrial matrix is labeled with a quickly folding, bright red fluorescent protein in a BacMam vector for controlled expression in a wide variety of cell types.	5 mL	\$595
#P1020R	CMV	Nuclear-targeted Montana Paintbrush (NLSPaint)	A quickly folding bright red fluorescent protein tagged with a nuclear localization signal labeling the nucleus in a BacMam vector for controlled expression in a wide variety of cell types.	5 mL	\$595
#P1020R	CMV	Nuclear-targeted Montana Paintbrush (NLSPaint)	A quickly folding bright red FP tagged with a nuclear localization signal labeling the nucleus in a BacMam vector for controlled expression in a wide variety of cell types.	5 mL	\$595
#P1030R	CMV	Lysosome-targeted Montana Paintbrush (Red Lamp2Paint)	Montana Paintbrush red FP targeted to the Lysosome with Lamp2. Packaged in our Big Sky BacMam vector for consistent, titratable transduction into a wide range of cell types	5 mL	\$595
#P1040R	CMV	Golgi-targeted Montana Paintbrush (Red GiantinPaint)	Montana Paintbrush red FP targeted to the Golgi with Giantin. Packaged in our Big Sky BacMam vector for consistent, titratable transduction into a wide range of cell types.	5 mL	\$595
#P1040G	CMV	Golgi-targeted mNeon Green (Green GiantinPaint)	mNeonGreen FP targeted to the Golgi with Giantin. Packaged in our Big Sky BacMam vector for consistent, titratable transduction into a wide range of cell types.	5 mL	\$595
#P1050R	CMV	Tau-targeted Montana Paintbrush (Red TauPaint)	Red TauPaint is the Montana Paintbrush red FP fused to tau. Packaged in our Big Sky BacMam vector for consistent, titratable transduction into a wide range of cell types.	5 mL	\$595
#P1050G	CMV	Tau-targeted mNeonGreen (Green TauPaint)	mNeonGreen FP fused to Tau. Packaged in our Big Sky BacMam vector for consistent, titratable transduction into a wide range of cell types.	5 mL	\$595
#P1060G	CMV	F-actin targeted mNeon Green (Green β -ActinPaint)	mNeonGreen FP fused to Human β -Actin. Packaged in our Big Sky BacMam vector for consistent, titratable transduction into a wide range of cell types.	5 mL	\$595
#P1060R	CMV	F-actin Targeted Montana Paintbrush (Red BetaActinPaint)	Montana Paintbrush red FP fused to human β -Actin. Packaged in our Big Sky BacMam vector for consistent, titratable transduction into a wide range of cell types	5 mL	\$595
#P1070G	CMV	Alpha Tubulin targeted mNeonGreen (Green AlphaTubulinPaint)	mNeonGreen FP fused to Alpha Tubulin. Packaged in our Big Sky BacMam vector for consistent, titratable transduction into a wide range of cell types.	5 mL	\$595
#P1070R	CMV	Alpha Tubulin targeted Montana Paintbrush (Red AlphaTubulinPaint)	Montana Paintbrush red FP fused to Alpha Tubulin. Packaged in our Big Sky BacMam vector for consistent, titratable transduction into a wide range of cell types.	5 mL	\$595
#P1080G	CMV	Endosome targeted mNeon Green (Green EndoPaint)	mNeonGreen FP targeted to the Endosome with FYVE. Packaged in our Big Sky BacMam vector for consistent, titratable transduction into a wide range of cell types.	5 mL	\$595
#P1090G	CMV	Omeasome-targeted mNeon Green (Green OmegaPaint)	mNeonGreen FP targeted to the Omeasome with ZNFN2A1. Packaged in our Big Sky BacMam vector for consistent, titratable transduction into a wide range of cell types.	5 mL	\$595

Product Code	Promoter	Neurodegenerative Associated Proteins	Use	Sizes	List Price (USD)
#N0100N #N0110N	CMV	Rhodopsin WT Rhodopsin P23H	Rhodopsin (Wild Type or P23H Mutation) in BacMam for expression in mammalian cells. The P23H mutation is associated with Retinitis Pigmentosa.	5 mL	\$395
#N0200N #N0210N	CMV	TDP-43 WT TDP-43 M337V	TDP-43 (Wild Type or M337V Mutation) in BacMam for expression in mammalian cells. The TDP-43 protein is associated with amyotrophic lateral sclerosis (ALS) and frontotemporal dementia.	5 mL	\$395
#N0300N #N0310N	CMV	FUS WT FUS P525L	FUS (Wild Type or P525L Mutation) packaged in BacMam for expression in mammalian cells. FUS is associated with amyotrophic lateral sclerosis (ALS) and frontotemporal dementia..	5 mL	\$395
#N0400N #N0410N #N0420N #N0430N #N0440N #N0450N	CMV	α -synuclein WT α -synuclein A53T α -synuclein A30P α -synuclein E46K α -synuclein H50Q α -synuclein G51D	α -Synuclein (Wild Type or Mutation) packaged in BacMam for expression in mammalian cells. α -Synuclein is associated with Parkinson's disease.	5 mL	\$395
#N0500N #N0510N #N0520N	CMV	SOD1 WT SOD1 G85R SOD1 D90A	SOD1 (Wild Type or Mutation) packaged in BacMam for expression in mammalian cells. SOD1 is associated with amyotrophic lateral sclerosis (ALS).	5 mL	\$395

Product Code	Promoter	GPCR Receptors in BacMam	Use	Sizes	List Price (USD)
#Z0100N	CMV	Human D ₁ Dopamine Receptor (DRD1)	Gs	30 mL	\$795
#Z0110N	CMV	Human D ₂ Dopamine Receptor (DRD2)	Gi	30 mL	\$795
#Z0120N	CMV	Human D ₃ Dopamine Receptor (DRD3)	Gi	30 mL	\$795
#Z0200N	CMV	Human M ₁ Muscarinic Receptor (CHRM1)	Gq	30 mL	\$795
#Z0200G	CMV	Human M ₁ Muscarinic Receptor (CHRM1) - Green	Gq	30 mL	\$795
#Z0200R	CMV	Human M ₁ Muscarinic Receptor (CHRM1) - Red	Gq	30 mL	\$795
#Z0210N	CMV	Human M ₂ Muscarinic Receptor (CHRM2)	Gi	30 mL	\$795
#Z0220N	CMV	Human M ₃ Muscarinic Receptor (CHRM3)	Gq	30 mL	\$795
#Z0300N	CMV	Human A ₁ Adenosine Receptor (ADORA1)	Gi	30 mL	\$795
#Z0400N	CMV	Human Succinate Receptor (SUCNR1)	Gi	30 mL	\$795
#Z0500N	CMV	Human β ₂ Adrenergic Receptor (ADRB2)	Gs	30 mL	\$795
#Z0500G	CMV	Human β ₂ Adrenergic Receptor (ADRB2) - Green	Gs	30 mL	\$795
#Z0500R	CMV	Human β ₂ Adrenergic Receptor (ADRB2) - Red	Gs	30 mL	\$795
#Z0600N	CMV	Human Glucagon-like Peptide GLP-1 Receptor (GLP1R)	Gs	30 mL	\$795
#Z0610N	CMV	Human Calcitonin Receptor (CALCR)	Gs/Gq	30 mL	\$795
#Z0620N	CMV	Human Vasopressin Receptor (AVPR2)	Gs	30 mL	\$795
#Z0700N	CMV	Human Nociceptin Opioid Receptor (OPRX)	Gi	30 mL	\$795
#Z0710N	CMV	Human Kappa Opioid Receptor (OPRK1)	Gi	30 mL	\$795
#Z0720N	CMV	Human Mu Opioid receptor (OPRM1)	Gi	30 mL	\$795
#Z0730N	CMV	Human Cannabinoid Receptor Type 1 (CNR1)	Gi	30 mL	\$795
#Z0740N	CMV	Human Delta Opioid Receptor (OPRD1)	Gi	30 mL	\$795
#Z0800N	CMV	Human Angiotensin II Type I Receptor (AGTR1)	Gq	30 mL	\$795
#Z0900N	CMV	Human Oxytocin Receptor (OXTR)	Gq	30 mL	\$795
#Z1000N	CMV	Human C-X-C Chemokine Receptor Type 4 (CXCR4)	Gi	30 mL	\$795
#Z1100N	CMV	Human Serotonin 5-HT _{2A} Receptor (HTR2A)	Gq	30 mL	\$795
#Z1200N	CMV	Human Proteinase-Activated Receptor 2 (PAR2-F2RL1)	Gq	30 mL	\$795
#Z1300N	CMV	Human Lysophosphatidic Acid Receptor 1 (LPAR1)	Gq, Gi, G ₁₂	30 mL	\$795
#Z1400N	CMV	Human Parathyroid Hormone 1 Receptor (PTH1R)	Gs	30 mL	\$795
#Z1500N	CMV	Human Serotonin 5-HT _{1A} Receptor (HTR1A)	Gi	30 mL	\$795
#Z1600N	CMV	Human Gastric Inhibitory Polypeptide Receptor (GIPR)	Gs	30 mL	\$795
#Z1700N	CMV	Human Cannabinoid Receptor Type 2 (CNR2)	Gi	30 mL	\$795
#Z1800N	CMV	Human Serotonin 5-HT _{2C} Receptor (HTR2C)	Gq	30 mL	\$795
#Z1900N	CMV	Human Serotonin 5-HT _{1B} Receptor (HTR1B)	Gi	30 mL	\$795
#Z2000N	CMV	Human Glucagon Receptor (GCGR)	Gs	30 mL	\$795
#Z2100N	CMV	Human Somatostatin Receptor Type 2 (SSTR2)	Gi	30 mL	\$795



Product Code	Promoter	Optogenetic Toolkit	Use	Sizes	List Price (USD)
#V0100N	CMV	bPAC	bPAC blue-light activated adenylyl cyclase which, when exposed to blue light produces cAMP without any additional proteins or cofactors.	10 mL	\$595
#V0101N	CMV	Kir 2.1	An inwardly rectifying potassium channel, can be used to hyperpolarize cell types such as HEK 293	10 mL	\$595
#V0102N	CMV	Na Channel	Bacterial, voltage-sensitive Na channel for creating artificial action potentials in fibroblasts .	10 mL	\$595
#V0103N	CMV	cAMP-gated channel	cAMP-gated ion channel . This is a mutant version of a cGMP-gated channel that responds to cAMP instead.	10 mL	\$595
#C0500N	CMV				
#C0505N	CAG	Cre Recombinase	Expresses Cre	10 mL	\$395
#C0506N	Synapsin				

Product Code	Promoter	Other BacMam Products	Use	Sizes	List Price (USD)
#X0000N	CMV	Empty BacMam	Expresses no protein, use as a negative control.	5 mL	\$295
#X0250X	CMV	Gs Mutant	A constitutively active Gs can be added during the transduction process to increase cAMP (can be used as add-on to Gi Kit #X0200G)	5 mL	\$295
#F0500G #F0505G	CMV CAG	mNeonGreen BacMam	Test expression efficiency in your cells.	5 mL	\$295
#F0520G	CMV	Nuclear-Targeted mNeonGreen BacMam	Expresses Nuclear-targeted mNeonGreen Fluorescent Protein	5 mL	\$295
#F0500R #F0505R	CMV CAG	Red Fluorescent Protein BacMam	Test expression efficiency in your cells.	5 mL	\$295
#D0700G	CMV	ArcLight Voltage	Monitor electrical activity in excitable cells with changes in green fluorescence.	30 mL	\$1695
#K0002N	CMV	GPCR Kinase-2 (GRK2)	GPCR Kinase-2 packaged in BacMam for expression in mammalian cells.	10 mL	\$695
#K0003N	CMV	GPCR Kinase-3 (GRK3)	GPCR Kinase-3 packaged in BacMam for expression in mammalian cells.	10 mL	\$695
#K0005N	CMV	GPCR Kinase-5 (GRK5)	GPCR Kinase-5 packaged in BacMam for expression in mammalian cells.	10 mL	\$695
#R0001N	CMV	RAMP1 BacMam	Receptor activity modifying protein 1 (RAMP1) packaged in Big Sky BacMam for expression in mammalian cells.	5 mL	\$595
#R0002N	CMV	RAMP2 BacMam	Receptor activity modifying protein 2 (RAMP2) packaged in Big Sky BacMam for expression in mammalian cells.	5 mL	\$595
#R0003N	CMV	RAMP3 BacMam	Receptor activity modifying protein 3 (RAMP3) packaged in Big Sky BacMam for expression in mammalian cells.	5 mL	\$595

