

Product Code	Promoter	BacMam Version	Product Name	Use	Sizes	List Price (USD)
Arrestin Recruitment Biosensors						
#D2000G	CMV	OG	Borealis – V2R	Kinetic arrestin signaling assay through activation of Vasopressin. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2010G	CMV	OG	Borealis – AT1R	Kinetic arrestin signaling assay through activation of Angiotensin. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2020G	CMV	OG	Borealis – GLP-1R	Kinetic arrestin signaling assay through activation of GLP-1R. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2030G	CMV	OG	Borealis – MOR	Kinetic arrestin signaling assay through activation of μ -Opioid receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2040G	CMV	OG	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.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2050G	CMV	OG	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.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2060G	CMV	OG	Borealis – OTR	Kinetic arrestin signaling assay through activation of Oxytocin receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2070G	CMV	OG	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.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2080G	CMV	OG	Borealis – KOR	Kinetic arrestin signaling assay through activation of κ -Opioid receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2090G	CMV	OG	Borealis – PAR2	Kinetic arrestin signaling assay through activation of the PAR2 receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2100G	CMV	OG	Borealis – DOR	Kinetic arrestin signaling assay through activation of the DOR receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2110G	CMV	OG	Borealis – LPA1R	Kinetic arrestin signaling assay through activation of the LPA1 receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2120G	CMV	OG	Borealis – PTH1R	Kinetic arrestin signaling assay through activation of the PTH1 receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2130G	CMV	OG	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.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2140G	CMV	OG	Borealis – GCGR	Kinetic arrestin signaling assay through activation of the Glucagon receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2150G	CMV	OG	Borealis – MC4R	Kinetic arrestin signaling assay through activation of the Melanocortin 4 receptor. Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2160G	CMV	OG	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.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2170G	CMV	OG	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.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2180G	CMV	OG	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.	10mL	\$895
					30mL	\$1795
					60mL	\$2995
#D2190G	CMV	OG	Borealis – ACKR3	Kinetic arrestin signaling assay through activation of the Apelin Receptor (APJ/APLNR). Monitor arrestin-3 signaling on fluorescence plate readers or imaging systems.	10mL	\$895
					30mL	\$1795
					60mL	\$2995



Product Code	Promoter	BacMam Version	Product Name	Use	Sizes	List Price (USD)
cADDIs cAMP Biosensors						
#D0250G	CMV Synapsin	Big Sky	Green Down cADDIs cAMP	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.	5mL	\$895
#D0251G					15mL	\$1795
					30mL	\$2995
#U0250G	CMV Synapsin	Big Sky	Green Up cADDIs cAMP	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.	5mL	\$895
#U0251G					15mL	\$1795
					30mL	\$2995
#U0250R	CMV	Big Sky	Red Up cADDIs cAMP	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.	5mL	\$895
					15mL	\$1795
					30mL	\$2995
#X0250G	CMV	Big Sky	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.	5mL	\$895
					15mL	\$1795
					30mL	\$2995
#D0201G	CMV	OG	Green Cilia-Targeted cADDIs cAMP	Monitor changes in cAMP in the primary cilia with high N.A. fluorescence microscopy.	10mL	\$995
#D0211G					10mL	\$995
#D0221G	CMV	OG	Green Nucleus-Targeted cADDIs cAMP	Monitor changes in cAMP in the cell nucleus with fluorescence imaging.	10mL	\$995
#U0221R					10mL	\$995
#D0231G	CMV	OG	Green Membrane-Targeted fMP cADDIs cAMP	Flag tagged MyrPalm cADDIs targeted to the plasma membrane.	10mL	\$995
#D0241G					10mL	\$995
#U0241R	CMV	OG	Red Membrane-Targeted fS15 cADDIs cAMP	Flag tagged S15 cADDIs targeted to the plasma membrane (red upward).	10mL	\$995
#U0261G					10mL	\$995
#U0261G	CMV	Big Sky	AKAP18a-Targeted cADDIs cAMP	Measure changes in signalsomes, with cADDIs sensor attached directly to different AKAP proteins.	5mL	\$895
#U0262G					5mL	\$895
#U0263G	CMV	Big Sky	AKAP1-Targeted cADDIs cAMP	Measure changes in signalsomes, with cADDIs sensor attached directly to different AKAP proteins.	5mL	\$895
#U0264G					5mL	\$895
#U0266G	CMV	Big Sky	AKAP12-Targeted cADDIs cAMP	Measure changes in signalsomes, with cADDIs sensor attached directly to different AKAP proteins.	5mL	\$895
#U0267G					5mL	\$895
#U0267G	CMV	Big Sky	smAKAP-Targeted cADDIs cAMP	Measure changes in signalsomes, with cADDIs sensor attached directly to different AKAP proteins.	5mL	\$895

Product Code	Promoter	BacMam Version	Product Name	Use	Sizes	List Price (USD)
DAG, PIP ₂ , Calcium, cGMP Biosensors						
#D0350G	CMV	Big Sky	Green Down DAG	Screen Gq-coupled GPCRs to monitor increases in DAG on fluorecence plate readers or imaging systems. Packaged in our new vector yielding significantly improved transduction efficiency.	5mL	\$895
					15mL	\$1795
					30mL	\$2995
#U0350G	CMV	Big Sky	Green Up DAG	Screen Gq-coupled GPCRs to monitor increases in DAG on fluorecence plate readers or imaging systems. Packaged in our new vector yielding significantly improved transduction efficiency.	5mL	\$895
					15mL	\$1795
					30mL	\$2995
#D0300R	CMV	Big Sky	Red Down DAG	Monitor Gq-mediated changes in diacylglycerol on fluorecence imaging systems.	10mL	\$895
					30mL	\$1795
#U0350R	CMV	Big Sky	Red Up DAG	Monitor Gq-mediated changes in diacylglycerol on fluorecence imaging systems	10mL	\$895
					30mL	\$1795
#D0331G	CMV	OG	Green Membrane-Targeted CAAX DAG	DAG sensor targeted (CAAX) to the plasma.	30mL	\$1795
#D0450G	CMV	Big Sky	Green PIP2 Sensor	Monitor changes in PIP ₂ on fluorecence imaging systems	10mL	\$895
#D0450R	CMV	Big Sky	Red PIP2 Sensor	Monitor changes in PIP ₂ on fluorecence imaging systems	10mL	\$895
#U0650G	CMV	Big Sky	Green GECO Ca ²⁺	Green Ca ²⁺ indicator packaged in our new vector yielding significantly improved transduction efficiency.	5mL	\$895
					15mL	\$1795
					30mL	\$2995
#U0650R	CMV	Big Sky	Red GECO Ca ²⁺	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.	5mL	\$895
					15mL	\$1795
					30mL	\$2995
#D0800G	CMV	OG	Green Down GENie cGMP	Detect changes in cGMP with a genetically encoded nucleotide indicator (GENie).	10mL	\$895
					30mL	\$1795
* CAG promoter kits also available, only in 30mL						

Product Code	Promoter	BacMam Version	Product Name	Use	Sizes	List Price (USD)
Cell Stress Biosensors						
#U0921G	CMV	Big Sky	IRE1-XBP1 Cell Stress	Bright fluorescent assay to detect response and reversal of chemical or genetic ER stress and UPR.	5mL	\$895
					15mL	\$1795
					30mL	\$2995

Product Code	Promoter	BacMam Version	SARS-CoV-2 Reagents	Use	Sizes	List Price (USD)
SARS-CoV-2 Reagents						
#C1100G	CMV	OG	mNeonGreen Fluorescent ACE2	Controlled expression of green fluorescent ACE2 for SARS-CoV-2 drug screening.	5mL	\$595
#C1100R	CMV	OG	Red Fluorescent ACE2	Controlled expression of red fluorescent ACE2 for SARS-CoV-2 drug screening.	5mL	\$595
#C1100N	CMV	OG	Unlabeled ACE2	Controlled expression of (unlabeled) ACE2 for SARS-CoV-2 drug screening.	5mL	\$595
#C1110G	CMV	Spike Pseudotyped	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.	5mL	\$595
#C1120G	CMV	Spike Pseudotyped	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.	5mL	\$595
#C1123G	CMV	Spike Pseudotyped	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.	5mL	\$595
#C1130N	CMV	OG	TMPRSS2	TMPRSS2 in BacMam for controlled expression in live mammalian cells and increased spike binding.	5mL	\$595
#C1140R	CMV	OG	Neuropilin-1 - Red	Red fluorescent labeled Neuropilin 1 in a BacMam vector for controlled expression in a wide variety of cell types	5mL	\$595
#C1140N	CMV	OG	Neuropilin-1 - Unlabeled	Unlabeled Neuropilin 1 in a BacMam vector for controlled expression in a wide variety of cell types.	5mL	\$595
#C1150N	CMV	OG	Cathepsin-L	Cathepsin L in BacMam for efficient delivery to practically any cell type.	5mL	\$595
#C1160G	CMV	OG	3CLglow	Green fluorescent biosensor for the SARS-CoV-2 Main protease. Screen for potential inhibitors of SARS-CoV-2 in living cells.	5mL	\$595
#C1161G	CMV	OG	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.	5mL	\$595
#C1162G	CMV	OG	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.	5mL	\$595
#C1200N	CMV	OG	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.	5mL	\$595
#C1201N	CMV	OG	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.	5mL	\$595
#C1202N	CMV	OG	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.	5mL	\$595
#C1203N	CMV	OG	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.	5mL	\$595
#C1204N	CMV	OG	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.	5mL	\$595
#C1205N	CMV	OG	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.	5mL	\$595
#C1206N	CMV	OG	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.	5mL	\$595
#C1207N	CMV	OG	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.	5mL	\$595
#C1208N	CMV	OG	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.	5mL	\$595
#C1209N	CMV	OG	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.	5mL	\$595
#C1210N	CMV	OG	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.	5mL	\$595
#C1211N	CMV	OG	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.	5mL	\$595
#C1212N	CMV	OG	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.	5mL	\$595
#C1213N	CMV	OG	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.	5mL	\$595
#C1214N	CMV	OG	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.	5mL	\$595
#C1215N	CMV	OG	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.	5mL	\$595



Product Code	Promoter	BacMam Version	Product Name	Use	Sizes	List Price (USD)
Cell Painting & Fluorescent Proteins						
#F0500G #F0505G	CMV CAG	OG	mNeonGreen BacMam	Test expression efficiency in your cells using our OG BacMam vector for expression in mammalian cells.	5mL	\$295
#P1000G #P1001G #P1005G #P1008G	CMV Synapsin CAG Ef1- α	Big Sky	mNeonGreen BacMam	Test expression efficiency in your cells using our Big Sky BacMam vector for consistent, titratable transduction into a wide range of cell types.	5mL	\$595
#P5000G	CMV	Evergreen	mNeonGreen BacMam	Test expression efficiency in your cells using our Evergreen BacMam vector for HDAC-inhibitor Free expression in human or primate cells..	5mL	\$595
#P1000R	CMV	Big Sky	Montana Paintbrush BacMam (MTPaint)	A quickly folding bright cytosolic red fluorescent protein.	5mL	\$595
#P1010G	CMV	Big Sky	Mitochondria-Targeted mNeonGreen (Green MitoPaint)	mNeonGreen FP targeted to the mitochondria with Cox8.	5mL	\$595
#P1010R	CMV	Big Sky	Mitochondrial-Targeted Montana Paintbrush (Red MitoPaint)	The inner mitochondrial matrix is labeled with a quickly folding, bright red fluorescent protein.	5mL	\$595
#P1020G	CMV	Big Sky	Nucleus-Targeted mNeonGreen (Green H2BPaint)	mNeonGreen FP targeted to the nucleus using H2B.	5mL	\$595
#P1020R	CMV	Big Sky	Nucleus-Targeted Montana Paintbrush (Red NLSPaint)	Montana Paintbrush RFP targeted to the nucleus of your cells, using NLS.	5mL	\$595
#P1030G	CMV	Big Sky	Lysosome-Targeted mNeonGreen (Green Lamp2Paint)	mNeonGreen fluorescent protein targeted to the Lysosome with Lamp2	5mL	\$595
#P1030R	CMV	Big Sky	Lysosome-Targeted Montana Paintbrush (Red Lamp2Paint)	Montana Paintbrush red FP targeted to the Lysosome with Lamp2.	5mL	\$595
#P1040G	CMV	Big Sky	Golgi-Targeted mNeonGreen (Green GiantinPaint)	mNeonGreen FP targeted to the Golgi with Giantin.	5mL	\$595
#P1040R	CMV	Big Sky	Golgi-Targeted Montana Paintbrush (Red GiantinPaint)	Montana Paintbrush red FP targeted to the Golgi with Giantin.	5mL	\$595
#P1050G	CMV	Big Sky	Tau-Targeted mNeonGreen (Green TauPaint)	mNeonGreen FP fused to Tau.	5mL	\$595
#P1050R	CMV	Big Sky	Tau-Targeted Montana Paintbrush (Red TauPaint)	Red TauPaint is the Montana Paintbrush red FP fused to tau..	5mL	\$595
#P1060G	CMV	Big Sky	F-actin-Targeted mNeonGreen (Green β -ActinPaint)	mNeonGreen FP fused to Human β -Actin.	5mL	\$595
#P1060R	CMV	Big Sky	F-actin-Targeted Montana Paintbrush (Red β -ActinPaint)	Montana Paintbrush red FP fused to human β -Actin.	5mL	\$595
#P1070G	CMV	Big Sky	Alpha Tubulin-Targeted mNeonGreen (Green AlphaTubulinPaint)	mNeonGreen FP fused to Alpha Tubulin..	5mL	\$595
#P1070R	CMV	Big Sky	Alpha Tubulin-Targeted Montana Paintbrush (Red AlphaTubulinPaint)	Montana Paintbrush red FP fused to Alpha Tubulin.	5mL	\$595
#P1080G	CMV	Big Sky	Endosome targeted mNeon Green (Green EndoPaint)	mNeonGreen FP targeted to the Endosome with FYVE.	5mL	\$595
#P1090G	CMV	Big Sky	Omegasome-targeted mNeon Green (Green OmegaPaint)	mNeonGreen FP targeted to the Omegasome with ZNFN2A1.	5mL	\$595

Product Code	Promoter	BacMam Version	Neurodegenerative Associated Proteins	Use	Sizes	List Price (USD)
Neurodegenerative Associated Proteins						
#N0100N #N0110N	CMV	OG	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.	30mL	\$995
#N0200N #N0210N	CMV	OG	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.	30mL	\$995
#N0300N #N0310N	CMV	OG	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..	30mL	\$995
#N0400N #N0410N #N0420N #N0430N #N0440N #N0450N	CMV	OG	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.	30mL	\$995
#N0500N #N0510N #N0520N	CMV	OG	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).	30mL	\$995

Product Code	Promoter	BacMam Version	Product Name	Use	Sizes	List Price (USD)
G-protein Coupled Receptors in BacMam						
#Z0100N	CMV	OG	Human D ₁ Dopamine Receptor (DRD1)	Gs	30mL	\$995
#Z0119N	CMV	Big Sky	Human D ₂ Dopamine Receptor (DRD2)	Gi	15mL	\$995
#Z0129N	CMV	Big Sky	Human D ₃ Dopamine Receptor (DRD3)	Gi	15mL	\$995
#Z0200N	CMV	OG	Human M ₁ Muscarinic Receptor (CHRM1)	Gq	30mL	\$995
#Z0200G	CMV	OG	Human M ₁ Muscarinic Receptor (CHRM1) - Green	Gq	30mL	\$995
#Z0200R	CMV	OG	Human M ₁ Muscarinic Receptor (CHRM1) - Red	Gq	30mL	\$995
#Z0219N	CMV	Big Sky	Human M ₂ Muscarinic Receptor (CHRM2)	Gi	15mL	\$995
#Z0220N	CMV	OG	Human M ₃ Muscarinic Receptor (CHRM3)	Gq	30mL	\$995
#Z0309N	CMV	Big Sky	Human A ₁ Adenosine Receptor (ADORA1)	Gi	15mL	\$995
#Z0400N	CMV	OG	Human Succinate Receptor (SUCNR1)	Gi	30mL	\$995
#Z0500N	CMV	OG	Human β2 Adrenergic Receptor (ADRB2)	Gs	30mL	\$995
#Z0500G	CMV	OG	Human β2 Adrenergic Receptor (ADRB2)- Green	Gs	30mL	\$995
#Z0500R	CMV	OG	Human β2 Adrenergic Receptor (ADRB2) - Red	Gs	30mL	\$995
#Z0609N	CMV	Big Sky	Human Glucagon-like Peptide GLP-1 Receptor (GLP1R)	Gs	15mL	\$995
#Z0610N	CMV	OG	Human Calcitonin Receptor (CALCR)	Gs/Gq	30mL	\$995
#Z0620N	CMV	OG	Human Vasopressin Receptor (AVPR2)	Gs	30mL	\$995
#Z0709N	CMV	Big Sky	Human Nociceptin Opioid Receptor (OPRX)	Gi	15mL	\$995
#Z0719N	CMV	Big Sky	Human Kappa Opioid Receptor (OPRK1)	Gi	15mL	\$995
#Z0729N	CMV	Big Sky	Human Mu Opioid receptor (OPRM1)	Gi	15mL	\$995
#Z0739N	CMV	Big Sky	Human Cannabinoid Receptor Type 1 (CNR1)	Gi	15mL	\$995
#Z0749N	CMV	Big Sky	Human Delta Opioid Receptor (OPRD1)	Gi	15mL	\$995
#Z0800N	CMV	OG	Human Angiotensin II Type I Receptor (AGTR1)	Gq	30mL	\$995
#Z0900N	CMV	OG	Human Oxytocin Receptor (OXTR)	Gq	30mL	\$995
#Z1000N	CMV	OG	Human C-X-C Chemokine Receptor Type 4 (CXCR4)	Gi	30mL	\$995
#Z1010N	CMV	OG	Human C-C Chemokine Receptor 5 (CCR5)	Gi	30mL	\$995
#Z1100N	CMV	OG	Human Serotonin 5-HT2A Receptor (HTR2A)	Gq	30mL	\$995
#Z1200N	CMV	OG	Human Protease-Activated Receptor 2 (PAR2-F2RL1)	Gq	30mL	\$995
#Z1219N	CMV	Big Sky	Human Protease-Activated Receptor 1 (PAR1)	Gq, G12, G13	15mL	\$995
#Z1300N	CMV	OG	Human Lysophosphatidic Acid Receptor 1 (LPAR1)	Gq, Gi, G12	30mL	\$995
#Z1400N	CMV	OG	Human Parathyroid Hormone 1 Receptor (PTH1R)	Gs	30mL	\$995
#Z1419N	CMV	Big Sky	Human Parathyroid Hormone 2 Receptor (PTH2R)	Gs	15mL	\$995
#Z1509N	CMV	Big Sky	Human Serotonin 5-HT1A Receptor (HTR1A)	Gi	15mL	\$995
#Z1600N	CMV	OG	Human Gastric Inhibitory Polypeptide Receptor (GIPR)	Gs	30mL	\$995
#Z1700N	CMV	OG	Human Cannabinoid Receptor Type 2 (CNR2)	Gi	30mL	\$995
#Z1800N	CMV	OG	Human Serotonin 5-HT2C Receptor (HTR2C)	Gq	30mL	\$995
#Z1909N	CMV	Big Sky	Human Serotonin 5-HT1B Receptor (HTR1B)	Gi	15mL	\$995
#Z2000N	CMV	OG	Human Glucagon Receptor (GCGR)	Gs	30mL	\$995
#Z2109N	CMV	Big Sky	Human Somatostatin Receptor Type 2 (SSTR2)	Gi	15mL	\$995
#Z2209N	CMV	Big Sky	Human Melanocortin 1 Receptor (MC1R)	Gs	15mL	\$995
#Z2229N	CMV	Big Sky	Human Melanocortin 3 Receptor (MC3R)	Gs	15mL	\$995
#Z2239N	CMV	Big Sky	Human Melanocortin 4 Receptor (MC4R)	Gs	15mL	\$995
#Z2309N	CMV	Big Sky	Human Apelin Receptor (APLNR)	Gi	15mL	\$995
#Z2409N	CMV	Big Sky	Human GPR17 Receptor	Gi	15mL	\$995
#Z2509N	CMV	Big Sky	Human Atypical Chemokine Receptor 3 (ACKR3)		15mL	\$995
#Z2609N	CMV	Big Sky	Calcitonin Receptor-like Receptor (CRLR or CALCRL)		15mL	\$995

Product Code	Promoter	BacMam Version	Product Name	Use	Sizes	List Price (USD)
Optogenetic Toolkit						
#V0100N	CMV	OG	bPAC	bPAC blue-light activated adenylyl cyclase which, when exposed to blue light produces cAMP without any additional proteins or cofactors.	30mL	\$995
#V0101N	CMV	OG	Kir 2.1	An inwardly rectifying potassium channel, can be used to hyperpolarize cell types such as HEK 293	30mL	\$995
#V0102N	CMV	OG	Na Channel	Bacterial, voltage-sensitive Na channel for creating artificial action potentials in fibroblasts.	30mL	\$995
#V0103N	CMV	OG	cAMP-gated channel	cAMP-gated ion channel. This is a mutant version of a cGMP-gated channel that responds to cAMP instead.	30mL	\$995
#C0500N	CMV	OG	Cre Recombinase	Expresses Cre	30mL	\$995
#C0505N	CAG					
#C0506N	Synapsin					

Product Code	Promoter	BacMam Version	Product Name	Use	Sizes	List Price (USD)
Other BacMam Products						
#X0000N	CMV	OG	Empty BacMam	Expresses no protein, use as a negative control.	5mL	\$395
#X0010N	CMV	Big Sky	Empty Big Sky BacMam	Big Sky BacMam that expresses no protein, use as a negative control.	5mL	\$395
#X0250X	CMV	OG	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)	30mL	\$995
#K0002N	CMV	OG	GPCR Kinase-2 (GRK2)	GPCR Kinase-2 packaged in BacMam for expression in mammalian cells.	10mL	\$695
#K0003N	CMV	OG	GPCR Kinase-3 (GRK3)	GPCR Kinase-3 packaged in BacMam for expression in mammalian cells.	10mL	\$695
#K0005N	CMV	OG	GPCR Kinase-5 (GRK5)	GPCR Kinase-5 packaged in BacMam for expression in mammalian cells.	10mL	\$695
#R0001N	CMV	Big Sky	RAMP1 BacMam	Receptor activity modifying protein 1 (RAMP1) packaged in Big Sky BacMam for expression in mammalian cells.	5mL	\$595
#R0002N	CMV	Big Sky	RAMP2 BacMam	Receptor activity modifying protein 2 (RAMP2) packaged in Big Sky BacMam for expression in mammalian cells.	5mL	\$595
#R0003N	CMV	Big Sky	RAMP3 BacMam	Receptor activity modifying protein 3 (RAMP3) packaged in Big Sky BacMam for expression in mammalian cells.	5mL	\$595