

Product Code	Promoter	Fluorescent Biosensor BacMam Kit	Use	Sizes	List Price (USD)
#C1100G	CMV	mNeonGreen Fluorescent ACE2	Controlled expression of green fluorescent ACE2 for SARS-CoV-2 drug screening.	5 mL	\$395
				30 mL	\$795
				60 mL	\$1195
				200 mL	A.Q.
#C1100R	CMV	Red Fluorescent ACE2	Controlled expression of red fluorescent ACE2 for SARS-CoV-2 drug screening.	5 mL	\$395
				30 mL	\$795
				60 mL	\$1195
				200 mL	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	\$595
				30 mL	\$2595
				60 mL	\$3895
				200 mL	A.Q.
#C1110R	CMV	Pseudo SARS CoV-2 Red Reporter	BacMam virus pseudotyped with SARS-CoV-2 spike protein that will produce red fluorescent nuclei upon successful entry into host cells expressing ACE2.	5 mL	\$595
				30 mL	\$2595
				60 mL	\$3895
				200 mL	A.Q.
#C1120G	CMV	Pseudo SARS CoV-2 D614G Mutation 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	\$595
				30 mL	\$2595
				60 mL	\$3895
				200 mL	A.Q.
#C1120R	CMV	Pseudo SARS CoV-2 D614G Mutation Red Reporter	BacMam virus pseudotyped with SARS-CoV-2 D614G mutated spike protein that binds with ACE2 and expresses red fluorescent protein at the nucleus 24 hours after viral entry	5 mL	\$595
				30 mL	\$2595
				60 mL	\$3895
				200 mL	A.Q.
#C1130N	CMV	TMPRSS2	TMPRSS2 in BacMam for controlled expression in live mammalian cells and increased spike binding	5 mL	\$395
				30 mL	\$795
				60 mL	\$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	\$395
				30 mL	\$795
#C1150N	CMV	Cathepsin-L	Cathepsin L in BacMam for efficient delivery to practically any cell type.	5 mL	\$395
				30 mL	\$795
#C1160G	CMV	3CLglow	The 3CLpro (Mpro) inhibitor assay kit produces bright green fluorescence in the presence of 3CL protease activity in living cells.	5 mL	\$595
				30 mL	\$2595
				60 mL	\$3895

Product Code	Promoter	Fluorescent Biosensor BacMam Kit	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	\$595
#D0205G	CAG			30 mL	\$1195
				60 mL	\$1795
				200 mL	A.Q.
#U0200G	CMV	Green Up cADDIs cAMP	Screen Gs-coupled GPCRs to monitor increases in cAMP on fluorescence plate readers or imaging systems	10 mL	\$595
#U0205G	CAG			30 mL	\$1195
				60 mL	\$1795
				200 mL	A.Q.
#U0200R	CMV	Red Up cADDIs cAMP	Screen Gs-coupled GPCRs to monitor increases in cAMP on fluorescence plate readers or imaging systems	10 mL	\$595
#U0205R	CAG			30 mL	\$1195
				60 mL	\$1795
#X0200G	CMV	Green Gi cADDIs cAMP	Screen Gi-coupled GPCRs to monitor cAMP reduction on fluorescence plate readers or imaging systems	10 mL	\$595
#X0205G	CAG			30 mL	\$1195
				60 mL	\$1795
				200 mL	A.Q.
#D0201G	CMV	Green Cilia-Targeted cADDIs cAMP	Monitor changes in cAMP in the primary cilia with high N.A. fluorescence microscopy.	10 mL	\$595
				30 mL	\$1195
#D0211G	CMV	Ratiometric Cilia-Targeted cADDIs cAMP	Monitor changes in cAMP in the primary cilia with high N.A. fluorescence microscopy.	10 mL	\$595
				30 mL	\$1195
#D0221G	CMV	Green Nucleus-Targeted cADDIs cAMP	Monitor changes in cAMP in the cell nucleus with fluorescence imaging.	10 mL	\$595
				30 mL	\$1195
#U0221R	CMV	Red Nucleus-Targeted cADDIs cAMP	Monitor changes in cAMP in the cell nucleus with fluorescence imaging.	10 mL	\$595
				30 mL	\$1195
#D0231G	CMV	Green Membrane-Targeted fMP cADDIs cAMP	Flag tagged MyrPalm cADDIs targeted to the plasma membrane.	10 mL	\$595
				30 mL	\$1195
#U0241R	CMV	Red Membrane-Targeted fS15 cADDIs cAMP	Flag tagged S15 cADDIs targeted to the plasma membrane. (red upward)	10 mL	\$595
				30 mL	\$1195
#D0241G	CMV	Green Membrane-Targeted fS15 cADDIs cAMP	Flag tagged S15 cADDIs targeted to the plasma membrane (green downward).	10 mL	\$595
				30 mL	\$1195
#D0300G	CMV	Green Down DAG	Screen Gq-coupled GPCRs to monitor increases in DAG on fluorescence plate readers or imaging systems	10 mL	\$595
#D0305G	CAG			30 mL	\$1195
				60 mL	\$1795
				200 mL	A.Q.
#U0300G	CMV	Green Up DAG	Screen Gq-coupled GPCRs to monitor increases in DAG on fluorescence plate readers or imaging systems	10 mL	\$595
#U0305G	CAG			30 mL	\$1195
				60 mL	\$1795
#D0308G	Ef1- α	Cre-Dependent Green Down DAG	Target expression to Cre-expressing cells to monitor diacylglycerol on fluorescence imaging systems	10 mL	\$595
				30 mL	\$1195
#D0300R	CMV	Red Down DAG	Monitor Gq-mediated changes in diacylglycerol on fluorescence imaging systems	10 mL	\$595
				30 mL	\$1195
				60 mL	\$1795
#U0300R	CMV	Red Up DAG	Monitor Gq-mediated changes in diacylglycerol on fluorescence imaging systems	10 mL	\$595
				30 mL	\$1195
#D0331G	CMV	Membrane-Targeted DAG	DAG sensor targeted (CAAX) to the plasma	10 mL	\$595
				30 mL	\$1195



Product Code	Promoter	Fluorescent Biosensor BacMam Kit	Use	Sizes	List Price (USD)
#D0400G #D0405G	CMV CAG	Green PIP2 Sensor	Monitor changes in PIP2 on fluorescence imaging systems	10 mL	\$595
#D0400R	CMV	Red PIP2 Sensor	Monitor changes in PIP2 on fluorescence imaging systems	10 mL	\$595
#U0600G #U0608G	CMV Ef1- α	Green GECO Ca ²⁺	Green Ca ²⁺ indicator with CMV or Ef1- α promoter.	10 mL 30 mL 60 mL 200 mL	\$595 \$1195 \$1795 A.Q.
#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 30 mL 60 mL 200 mL	\$595 \$1195 \$1795 A.Q.
#D0700G	CMV	ArcLight Voltage	Monitor electrical activity in excitable cells with changes in green fluorescence.	10 mL	\$595
#D0800G	CMV	Green Down GENie cGMP	Detect changes in cGMP with a genetically encoded nucleotide indicator (GENie).	10 mL	\$595
#U0900G	CMV	Green Cell Stress	Bright fluorescent assay to detect response and reversal of chemical or genetic ER stress and UPR.	10 mL 30 mL 60 mL 200 mL	\$595 \$1195
#U0901G #U0915G	CMV CAG	Green/Red Ratiometric Cell Stress	Bright fluorescent assay to detect response and reversal of chemical or genetic ER stress and UPR.	10 mL 30 mL 60 mL 200 mL	\$595 \$1195 \$1795 A.Q.
#D1000G	CMV	β -Arrestin – Vasopressin	Kinetic β -arrestin signaling assay through activation of Vasopressin. Monitor β -arrestin signaling on fluorescence plate readers or imaging systems.	10 mL 30 mL 60 mL	\$895 \$1795 \$2695
#D1010G	CMV	β -Arrestin – Angiotensin	Kinetic β -arrestin signaling assay through activation of Angiotensin. Monitor β -arrestin signaling on fluorescence plate readers or imaging systems.	10 mL 30 mL 60 mL	\$895 \$1795 \$2695
#D1020G	CMV	β -Arrestin – GLP 1R	Kinetic β -arrestin signaling assay through activation of GLP-1R. Monitor β -arrestin signaling on fluorescence plate readers or imaging systems.	10 mL 30 mL 60 mL	\$895 \$1795 \$2695
#D1030G	CMV	β -Arrestin – Mu Opioid	Kinetic β -arrestin signaling assay through activation of mu opioid receptor. Monitor β -arrestin signaling on fluorescence plate readers or imaging systems.	10 mL 30 mL 60 mL	\$895 \$1795 \$2695
#D1040G	CMV	β -Arrestin – β 2 Adrenergic	Kinetic β -arrestin signaling assay through activation of β 2-Adrenergic receptor. Monitor β -arrestin signaling on fluorescence plate readers or imaging systems.	10 mL 30 mL 60 mL	\$895 \$1795 \$2695

Product Code	Promoter	GPCR Receptors in BacMam	Use	Sizes	List Price (USD)
#Z0100N	CMV	D1 Dopamine Receptor	Gs	10 mL 30 mL	\$395 \$795
#Z0110N	CMV	D2 Dopamine Receptor	Gi	10 mL 30 mL	\$395 \$795
#Z0115N	CAG				
#Z0120N	CMV	D3 Dopamine Receptor	Gi	10 mL 30 mL	\$395 \$795
#Z0200N	CMV	Human M1 Muscarinic Receptor	Gq	10 mL 30 mL	\$395 \$795
#Z0205N	CAG				
#Z0200G	CMV	Human M1 Muscarinic Receptor - Green	Gq	10 mL 30 mL	\$395 \$795
#Z0200R	CMV	Human M1 Muscarinic Receptor - Red	Gq	10 mL 30 mL	\$395 \$795
#Z0210N	CMV	Human M2 Muscarinic Receptor	Gi	10 mL 30 mL	\$395 \$795
#Z0215N	CAG				
#Z0220N	CMV	Human M3 Muscarinic Receptor		10 mL 30 mL	\$395 \$795
#Z0300N	CMV	A1 Adenosine Receptor	Gi	10 mL 30 mL	\$395 \$795
#Z0400N	CMV	Succinate Receptor	Gi	10 mL 30 mL	\$395 \$795
#Z0500N	CMV	β 2 Adrenergic Receptor	Gs	10 mL 30 mL	\$395 \$795
#Z0505N	CAG				
#Z0500G	CMV	β 2 Adrenergic Receptor - Green	Gs	10 mL 30 mL	\$395 \$795
#Z0500R	CMV	β 2 Adrenergic Receptor - Red	Gs	10 mL 30 mL	\$395 \$795
#Z0600N	CMV	Human GLP-1 Receptor		10 mL 30 mL	\$395 \$795
#Z0610N	CMV	Calcitonin Receptor	Gs/Gq	10 mL 30 mL	\$395 \$795
#Z0620N	CMV	Human Vasopressin Receptor		10 mL 30 mL	\$395 \$795
#Z0700N	CMV	Human Nociceptin Opioid Receptor (NOP) (Gi)	Gs	10 mL 30 mL	\$395 \$795
#Z0710N	CMV	Human Kappa Opioid Receptor (OPRK1) (Gi)	Gi	10 mL 30 mL	\$395 \$795
#Z0720N	CMV	Mu Opioid receptor (OPRM1) (Gi)	Gi	10 mL 30 mL	\$395 \$795
#Z0730N	CMV	CB1 Cannabinoid Receptor (CNR1)	Gi	10 mL 30 mL	\$395 \$795
#Z0800N	CMV	Human Angiotensin II Type I Receptor		10 mL 30 mL	\$395 \$795



Product Code	Promoter	Neurodegenerative Associated Proteins	Use	Sizes	List Price (USD)
#N0100N #N0110N	CMV	Rhodopsin WT Rhodopsin P23H	Retinitis Pigmentosa	5 mL	\$295
#N0200N #N0210N	CMV	TDP-43 WT TDP-43 M337V	Amyotrophic lateral sclerosis / Frontotemporal dementia	5 mL	\$295
#N0300N #N0310N	CMV	FUS WT FUS P525L	Amyotrophic lateral sclerosis / Frontotemporal dementia	5 mL	\$295
#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	Parkinson's Disease	5 mL	\$295
#N0500N #N0510N #N0520N	CMV	SOD1 WT SOD1 G85R SOD1 D90A	Amyotrophic lateral sclerosis	5 mL	\$295

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 #C0505N #C0506N	CMV CAG Synapsin	Cre Recombinase	Expresses Cre	10 mL	\$395

Product Code	Promoter	Other BacMam Products	Use	Sizes	List Price (USD)
#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	mNeon Green BacMam	Test expression efficiency in your cells.	5 mL	\$295.
#F0500R	CMV	Red Fluorescent Protein BacMam	Test expression efficiency in your cells.	5 mL	\$295