

Product Code	Product Description	Sizes	~Assays in 96 Well	~Assays in 384 Well	List Price (USD)	Vector	Plate Reader	Availability
#D0200G	Green Downward cADDIs cAMP (CMV Promoter)	5 mL 10mL	200 400	660 1350	\$395 \$595	BacMam	Z' > 0.9	immediate
#D0205G	Green Downward cADDIs cAMP (CAG Promoter)	30 mL 60 mL	1200 2400	4000 8000	\$1195 \$1795			
#U0200G	Green Upward cADDIs cAMP (CMV Promoter)	5 mL 10mL	200 400	660 1350	\$395 \$595			
#U0205G	Green Upward cADDIs cAMP (CAG Promoter)	30 mL 60 mL	1200 2400	4000 8000	\$1195 \$1795			
#U0200R	Red Upward cADDIs cAMP (CMV Promoter)	5 mL 10mL	200 400	660 1350	\$395 \$595	BacMam	Z' > 0.8	immediate
#U0205R	Red Upward cADDIs cAMP (CAG Promoter)	30 mL 60 mL	1200 2400	4000 8000	\$1195 \$1795			
#X0200G	Green Gi cADDIs cAMP (CMV Promoter)	5 mL 10mL	200 400	660 1350	\$395 \$595			
#X0205G	Green Gi cADDIs cAMP (CAG Promoter)	30 mL 60 mL	1200 2400	4000 8000	\$1195 \$1795			
#D0201G	Green Cilia-Targeted cADDIs cAMP	5 mL 10mL 30 mL 60 mL	200 400 1200 2400	660 1350 4000 8000	\$395 \$595 \$1195 \$1795	BacMam	N/A	immediate
#D0211G	Ratiometric Cilia-Targeted Green cADDIs cAMP	5 mL 10mL 30 mL 60 mL	200 400 1200 2400	660 1350 4000 8000	\$395 \$595 \$1195 \$1795	BacMam	N/A	immediate
#D0300G	Green Downward DAG (CMV Promoter)	5 mL 10mL	200 400	660 1350	\$395 \$595	BacMam	Z' > 0.8	immediate
#D0305G	Green Downward DAG (CAG Promoter)	30 mL 60 mL	1200 2400	4000 8000	\$1195 \$1795			
#U0300G	Green Upward DAG (CMV Promoter)	5 mL 10mL	200 400	660 1350	\$395 \$595			
#U0305G	Green Upward DAG (CAG Promoter)	30 mL 60 mL	1200 2400	4000 8000	\$1195 \$1795			
#D0308G	Cre-Dependent Green Downward DAG (Ef1- $\alpha$ Promoter)	5mL 10mL 30 mL	200 400 1200	660 1350 4000	\$395 \$595 \$1195	BacMam	N/A	immediate
#D0300R	Red Downward DAG	5 mL 10mL 30 mL 60 mL	200 400 1200 2400	660 1350 4000 8000	\$395 \$595 \$1195 \$1795	BacMam	Z' > 0.7	immediate
#U0300R	Red Upward DAG	10mL 30 mL 60 mL	400 1200 2400	1350 4000 8000	\$595 \$1195 \$1795	BacMam	Z' > 0.65	immediate
#D0400G	Green PIP <sub>2</sub> Sensor (CMV Promoter)	10 mL	400	1350	\$595	BacMam	N/A	immediate
#D0405G	Green PIP <sub>2</sub> Sensor (CAG Promoter)							(#D0405G coming soon)
#D0400R	Red PIP <sub>2</sub> Sensor	10 mL	400	1350	\$595	BacMam	N/A	immediate

Special promoter systems and targeted expression to organelles are [available by request](#).

[Ask us about](#) Lenti and AAV vectors or Cre-dependent expression

**List Price for US & Canada, Effective: January 1, 2017**



Product Code	Product Description	Sizes	~Assays in 96 Well	~Assays in 384 Well	List Price (USD)	Vector	Plate Reader	Availability
#C0500N	Cre Recombinase (CMV Promoter)	10 mL	-	-	\$495	BacMam	N/A	immediate
#C0505N	Cre Recombinase (CAG Promoter)							
#U0600R	Red GECO Ca <sup>2+</sup> Sensor	5 mL 10mL 30 mL 60 mL	200 400 1200 2400	660 1350 4000 8000	\$395 \$595 \$1195 \$1795	BacMam	Z' = 0.5	immediate
#D0700G	ArcLight Voltage Sensor	10 mL	400	1350	\$595	BacMam	N/A	immediate
#D0701G	Synapsin promoter ArcLight Voltage Sensor	10 mL	400	1350	\$595	BacMam	N/A	immediate
#D0800G	Green Downward cGMP	10 mL	400	1350	\$595	BacMam	N/A	early state Aug 1
#U0900G	Green Cell Stress Sensor	10mL 30mL 60mL	400 1200 2400	1350 4000 8000	\$595 \$1195 \$1795	BacMam	N/A	immediate
#U0901G	Red NLS + 2A + Green Stress Sensor	10mL 30mL 60mL	400 1200 2400	1350 4000 8000	\$595 \$1195 \$1795	BacMam	N/A	immediate