



6042 Cornerstone Court West, Suite B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### **FAM-cGMP Substrate Catalog #: 60201**

**DESCRIPTION:** The fluorogenic PDE substrate 2 {2'-(6-[Fluoresceineinyl] aminoethylcarbamoyl)guanosine-3',5'-cyclic monophosphate} is used for phosphodiesterase assays. It has been used in activity assays for PDE5, PDE6 and PDE9.

**PURITY:** >95% by HPLC.

**APPLICATIONS:** Study for phosphodiesterase activity, and screening of small molecule inhibitors for drug discovery and HTS applications.

**SUPPLIED AS:** 1 mM Aqueous solution

**QUANTITY:** 100  $\mu$ l

**STORAGE:** -20°C or -70°C.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)