



11526 Sorrento Valley Rd SteA2  
San Diego Ca 92121  
Tel: 1.858.202.1401  
1.858.829.3082  
Fax: 1.858.259.5734  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

***PDE3A*** (human, recombinant, N-terminal GST tag)  
Catalog #: 60030

**Description:** Human PDE3A (GenBank Accession No. NM\_000921), (a.a. 484-1141) with N-terminal GST tag, MW= 103.3 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

**Source:** Baculovirus infected Sf9 cells

**Formulated in:** 25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20, 50% glycerol, 20 mM glutathione, and 3 mM DTT.

**Stability:** >6 months at -80°C

**Purity:** >50% by SDS-PAGE.

**Specific Activity:** 1380 U/ug. 1 unit is defined as the amount of enzyme that will convert 1 pmole of 3', 5'-cAMP to 5' AMP per min at 30 °C in a reaction buffer. Assay buffer: 10 mM Tris-HCl, pH7.4, 10 mM MgCl<sub>2</sub>, 1 mM MnCl<sub>2</sub>, 0.4 mM 3', 5'-cAMP substrate, and 1 ng/μl PDE3A.

**Application:** useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### **Reference:**

1. Hung,S.H., Zhang,W., Pixley,R.A., Jameson,B.A., Huang,Y.C., Colman,R.F. and Colman,R.W. J. Biol. Chem. 281 (39), 29236-29244 (2006)
2. Hambleton,R., Krall,J., Tikishvili,E., Honeggar,M., Ahmad,F., Manganiello,V.C. and Movsesian,M.A. J. Biol. Chem. 280 (47), 39168-39174 (2005)

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** or **1.858.829.3082**, Fax **1.858.259.5734**  
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)