

## Data Sheet

### **PDE5A Mouse**

Mouse, recombinant, N-terminal GST tag

**Catalog #:** 60051

Lot #: 80207

Conc.: 0.34 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0,  
 124 mM NaCl, 2.4 mM KCl, 3 mM DTT, 18  
 mM glutathione, 10% glycerol.

**Stability:** >6 months at -80°C

**References:**

1. Das, A. *et al.*, *J. Biol. Chem.* **280** (13), 12944-12955 (2005).
2. Champion, H.C. *et al.*, *Proc. Natl. Acad. Sci. U.S.A.* **102** (5), 1661-1666 (2005).

**Description:**

Mouse PDE5A (GenBank Accession No. NM\_153422), full length, with N-terminal GST tag, MW=124 kDa, expressed in a Baculovirus-infected Sf9 cell expression system.

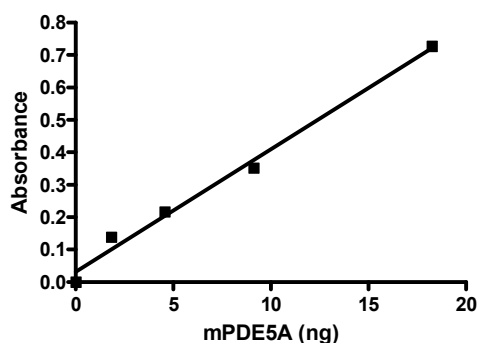
**Specific Activity:** 4630 pmol/min/μg.  
 Assay conditions: 10 mM Tris-HCl, pH 7.4,  
 10 mM MgCl<sub>2</sub>, 1 mM MnCl<sub>2</sub>, 0.1 mg/ml  
 BSA, 200 μM cGMP, 2.5 kU 5' nucleotidase,  
 37°C, 20 min.

**Application:**

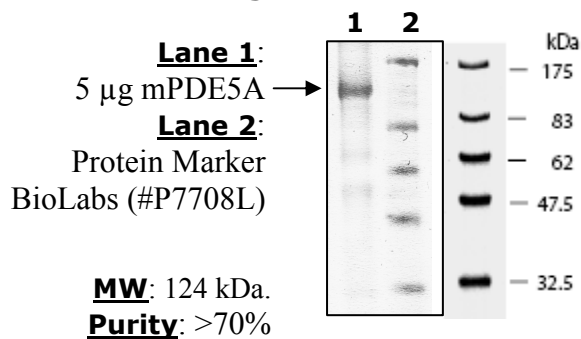
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### **Quality Assurance**

**mPDE5A Specific Activity**



**10% SDS-PAGE  
 Coomassie staining**



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

To place your order, please contact us by Phone **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)