

## Data Sheet

### PDE2A V1 (Mouse), GST-tag

Mouse, Recombinant, N-terminal GST-tag

**Catalog #:** 60017

**Lot #:** 160902-G1      **Conc.:** 0.21 mg/ml

**Formulated in:** 40mM Tris-HCl, pH 8.0, 110mM NaCl, 2.2mM KCl, 0.04% Tween-20 and 20% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

- Sadhu, K., *et al.*, *J. Histochem. Cytochem.* **47(7)**, 889-894 (1999).
- Iffland, A., *et al.*, *Biochemistry* **44 (23)**, 8312-8325 (2005).

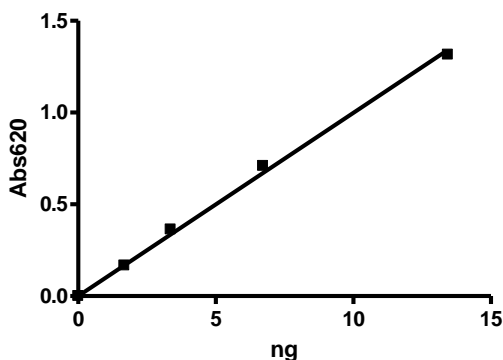
**Description:** Murine PDE2A, variant 1 (GenBank Accession No. NM\_001143848), a.a 2-939 (end), with N-terminal GST tag, MW=132 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

**Specific Activity:**  $\geq 20,000$  pmole/min/ $\mu\text{g}$ . One unit is defined as the amount of enzyme that will convert 1 pmole of 3', 5'-cAMP to 5'-AMP per min at  $37^{\circ}\text{C}$ . Assay condition: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl<sub>2</sub>, 0.1 mg/ml BSA, 0.05% Tween-20, 200  $\mu\text{M}$  cAMP, and serial dilutions of mouse PDE2A, mouse PDE4C and mouse PDE1B at  $37^{\circ}\text{C}$  for 20 min. Quantified by 5' nucleotidase cleaving the 5'-AMP product and releasing the phosphate group which is detected by Malachite Green Reagent.

**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity

## Quality Assurance

**Specific Activity**

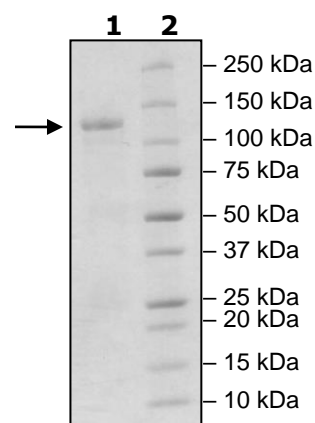


**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
2 $\mu\text{g}$  mPDE2A(v1)

**Lane 2:**  
Protein Marker

**MW:** 132 kDa  
**Purity:**  $\geq 86\%$



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 **1.858.481.8694**  
 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)