

Data Sheet

PDE11A, GST-Tag

Mouse, Recombinant, N-terminal GST tag

Catalog #: 60064

Lot #: 161205-1 **Conc.:** 0.4 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 20 mM Glutathione, 10% glycerol, and 3 mM DTT.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

References:

1. Fawcett, L. *et al.*, *PNAS*. **97(7)**, 3207-3707 (2000).
2. Horvath, A. *et al.*, *Nature Genetics*. **38**, 794-800 (2006).

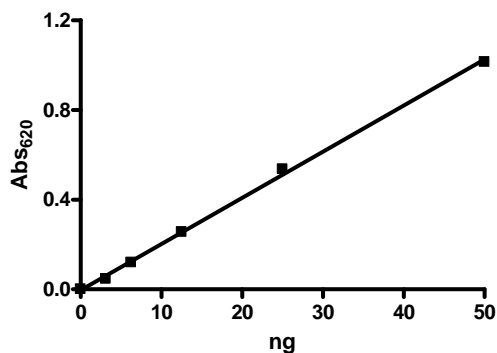
Description: Mouse PDE11A (GenBank Accession No. NM_001081033), a.a. 2-933(end) with N-terminal GST tag, MW= 131 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: $\geq 1560 \text{ pmol/min}/\mu\text{g}$. 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°C . Assay condition: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl_2 , 0.1 mg/ml BSA, 0.05% Tween-20, 200 μM cAMP, and serial dilutions of PDE11A at 37°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 profiling.

Quality Assurance

Specific Activity



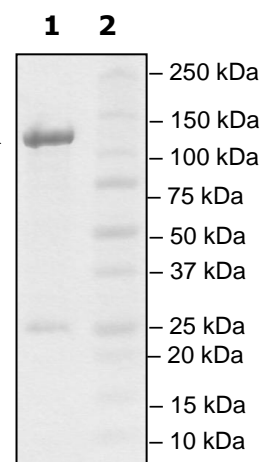
4-20% SDS-PAGE Coomassie staining

Lane 1:
PDE11A →

Lane 2:
Protein Marker

Purity: $\geq 85\%$

MW: 131 kDa



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

Please visit our website at: www.bpsbioscience.com