

Data Sheet

PDE9A, isoform b, GST-tag

Human, Recombinant, N-terminal GST tag
Catalog #: 60090
Lot #: 150831-G **Conc.:** 0.36 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 3 mM DTT,
 0.04% Tween-20 and 20% Glycerol.

Stability: At least 6 months at -80°C .
 Avoid freeze/thaw cycles. Storing diluted
 enzyme is not recommended, if
 necessary, use carrier protein (BSA 0.1 –
 0.5%).

References:

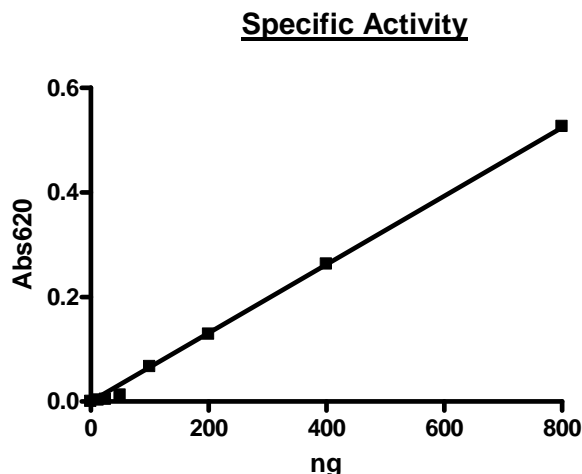
1. Rentero C. and Puigdomenech P. *BMC Mol. Biol.* **7**, 39 (2006).
2. Huai Q. *et al.*, *Proc. Natl. Acad. Sci. U.S.A.* **101** (26), 9624-9629 (2004).

Description: Human PDE9A, isoform b, also known as PDE9A2 (GenBank Accession No. NM_001001567), full length with N-terminal GST tag, MW=91 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: ≥ 200 pmol/min/ μg . Assay conditions: One unit is defined as the amount of enzyme that will convert 1 pmole of 3', 5'-cGMP to 5'-GMP per min at 37°C . Assay condition: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl₂, 0.1 mg/ml BSA, 0.05% Tween-20, 200 μM cGMP, and serial dilutions of PDE9A2 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.

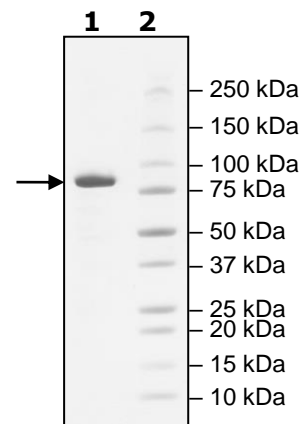
Quality Assurance



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

Lane 1:
 2 μg PDE9A2
Lane 2:
 Protein Marker

MW: 91 kDa
Purity: $\geq 92\%$



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