

Data Sheet

PDE7B

Rat, recombinant, N-terminal GST-tag

Catalog #: 60075

Lot #: 101117 **Conc.:** 0.25 mg/ml

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

Stability: >6 months at -80°C

References:

1. Gardner, C., *et al.* (2000). *Biochem Biophys Res Commun.* May 27;**272(1)**:186-92.
2. Reyes-Irisarri, E., *et al.* (2005). *Neuroscience* **132(4)**:1173-85.
3. Zhang, L., *et al.* (2008). *Proc. Natl. Acad. Sci. USA* Dec 9;**105(49)**:19532-7.

Description:

Rat PDE7B (GenBank Accession No. NM_080894), a.a. 108-446(end), with N-terminal GST-tag, MW=65 kDa, expressed in a Baculovirus-infected Sf9 cell expression system.

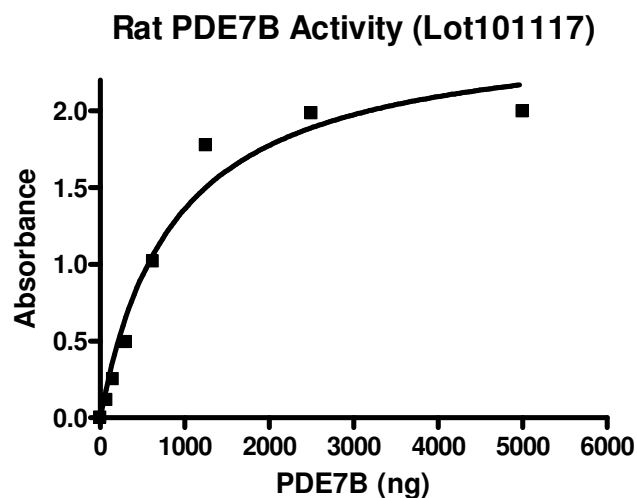
Specific Activity: 324 pmol/min/μg.

Assay condition: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl₂, 0.05% Tween-20, 200 μM cAMP, 0.1 mg/ml BSA, and 0.5-16 ng/μl rat PDE7B 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



10% SDS-PAGE Coomassie staining

Lane 1:

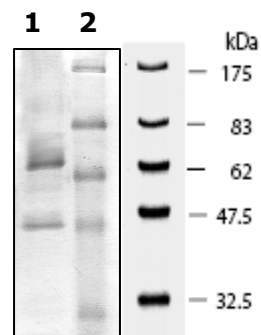
4 μg PDE7B →

Lane 2:

Protein Marker
 BioLabs (#P7708L)

MW: 65 kDa.

Purity: 30%



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