



11526 Sorrento Valley Rd SteA2
San Diego Ca 92121
Tel: 1.858.202.1401
1.858.829.3082
Fax: 1.858.259.5734
Email: info@bpsbioscience.com

Data Sheet

PDE7B

human, recombinant, , N-terminal GST tag
Catalog #: 60071

Formulated in: 25 mM Tris-HCl, pH 8.0,
100 mM NaCl, 0.05% Tween-20, 50% glycerol,
and 3 mM DTT.

Stability: >6 months at -80°C

References:

1. Gardner,C., et al., Biochem. Biophys. Res. Commun. 272 (1), 186-192 (2000)
2. Sasaki,T., et al., Biochem. Biophys. Res. Commun. 271 (3), 575-583 (2000)

Description:

Human PDE7B GenBank Accession No. NM_018945), (a.a. 107-end) with N-terminal GST tag, MW=66 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

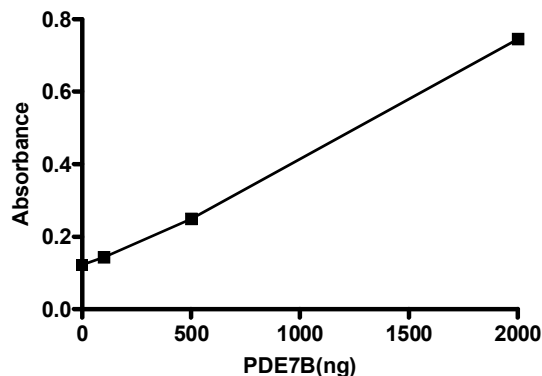
Specific Activity: 53 pmol/min/μg. 1 unit is defined as the amount of enzyme that will convert 1 pmole of 3', 5'-cAMP to 5' AMP per min at 30 °C in a reaction buffer. Assay conditions: 10 mM Tris-HCl, pH7.4, 10 mM MgCl₂, 1 mM MnCl₂, 400 μM cAMP, 2.5 kU 5' nucleotidase, 37°C, 20 min.

Application:

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

Quality Assurance

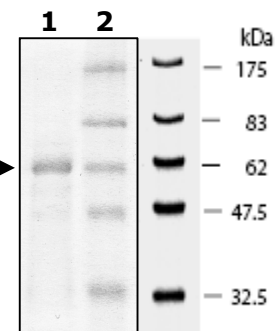
PDE7B Specific Activity



10% SDS-PAGE Coomassie staining

Lane 1:
4 μg PDE7B →

Lane 2:
Protein Marker
BioLabs (#P7708L)



MW: 66 kDa.
Purity: >80%

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

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com