



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

PDE7A1

human, recombinant, , N-terminal GST tag
Catalog #: 60070

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. Wang,H., et al., J. Biol. Chem. 280 (35), 30949-30955 (2005)
2. Torras-Llort,M. and Azorin,F., Biochem. J. 373 (PT 3), 835-843 (2003)

Description:

Human PDE7A1 GenBank Accession No. NM_002603), (a.a. 121-end) with N-terminal GST tag, MW=66 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 350 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.

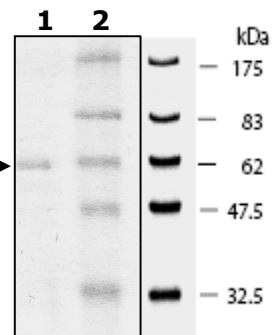
Quality Assurance

**10% SDS-PAGE
Coomassie staining**

Lane 1:
1 μg PDE7A

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