

Data Sheet

PDE3A (669-end)

Human, recombinant, N-terminal His-GST tag
Catalog #: 60032
Lot#: 110802HG **Conc.:** 0.46 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 16 mM
 glutathione, 20% Glycerol and 3 mM DTT.

Stability: >6 months at -80°C

References:

1. Hung, S.H., et al. *J. Biol. Chem.* **281** (39), 29236-29244 (2006)
2. Hambleton, R., et al. *J. Biol. Chem.* **280** (47), 39168-39174 (2005)

Description:

Human PDE3A (GenBank Accession No. NM_000921), amino acids 669-1141(end), with N-terminal His-GST tag, MW=82 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

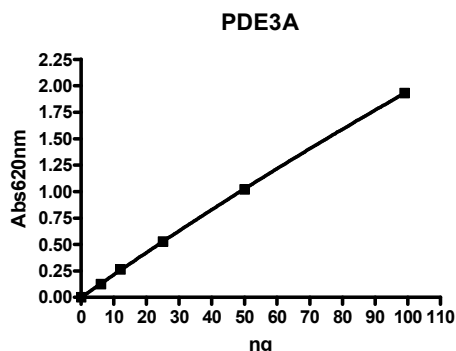
Specific Activity: 4000 pmol/min/μg.

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: 200 μM cAMP for 20 minutes at 37°C, followed by addition of Malachite Green reagents for 10 minutes at room temperature.

Application:

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

Quality Assurance

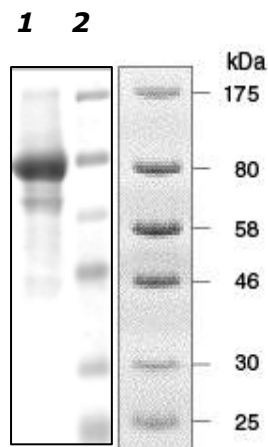


**10% SDS-PAGE
 Coomassie staining**

Lane 1:
 5 μg PDE3A

Lane 2:
 Protein Marker
 BioLabs (#P7708L)

MW: 82 kDa
Purity: ≥82%



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