

Data Sheet

PDE3A (Mouse), GST-Tag

Mouse, Recombinant, N-terminal GST-tag
Catalog #: 60036
Lot #: 170119-3 **Conc.:** 0.43 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol, 18 mM Glutathione

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:

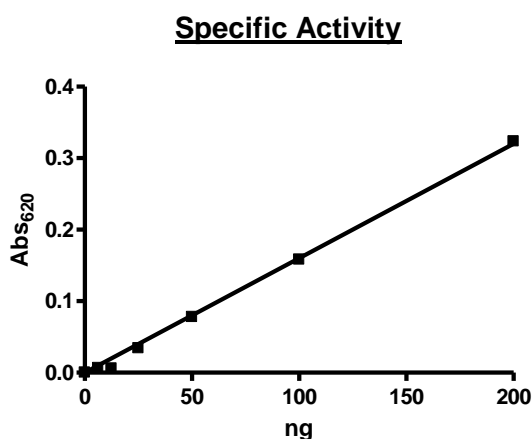
1. Hung, S.H., *et al.* *J. Biol. Chem.* **281** (39), 29236-29244 (2006)
2. Hambleton, R., *et al.* *J. Biol. Chem.* **280** (47), 39168-39174 (2005)

Description: Mouse PDE3A, GenBank Accession No. NM_018779, amino acids 484-1141 (end), with N-terminal GST tag, MW = 100 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: >280 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: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl₂, 0.1 mg/ml BSA, 0.05% Tween-20, 200 μM cAMP, and serial dilutions of Mouse PDE3A 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

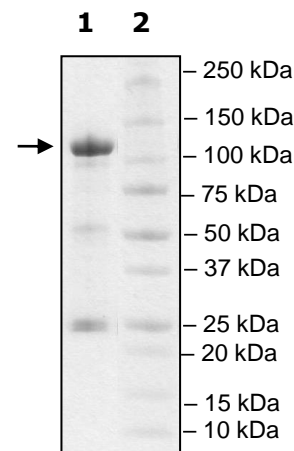


4-20% SDS-PAGE
Coomassie staining

Lane 1:
PDE3A(484-e)

Lane 2:
Protein Marker

MW: 100 kDa
Purity: $\geq 70\%$



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