

## Data Sheet

### PDE3B

Human, recombinant, N-terminal GST tag

**Catalog #:** 60031

**Lot#:** 110215G2      **Conc.:** 0.067 mg/ml

**Formulated in:** 40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 0.04% Tween-20, 20% glycerol

**Stability:** >6 months at -80 °C

### References:

1. Raymond, D.R. *et al.*, *Cell. Signal.* **19** (12), 2507-2518 (2007).
2. Palmer, D. *et al.*, *J. Biol. Chem.* **282** (13), 9411-9419 (2007).

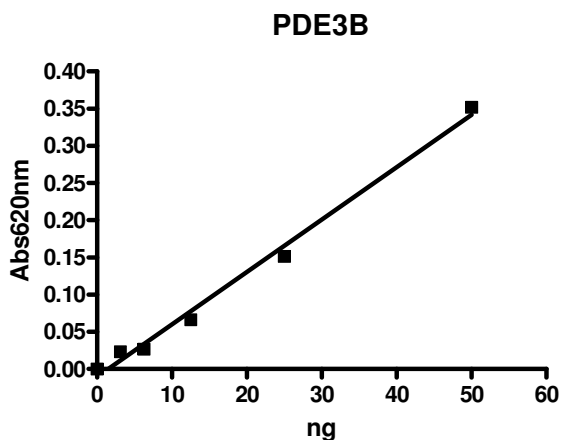
**Description:** Human PDE3B (GenBank Accession No. NM\_000922), (amino acids 592-end), with N-terminal GST tag, MW = 86 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

**Specific Activity:** 1270 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 MgCl<sub>2</sub>, 1 mM MnCl<sub>2</sub>, 200 μM cAMP, and 0.2 ng/μl PDE3B 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

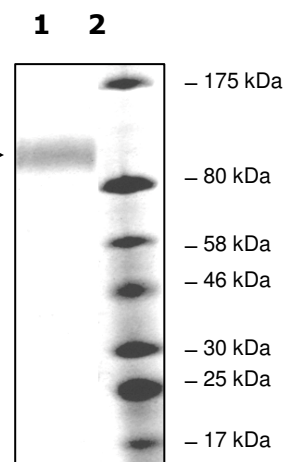


### 10% SDS-PAGE Coomassie staining

**Lane 1:**  
4 μg PDE3B

**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 86 kDa  
**Purity:** ≥90%



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)