

## Data Sheet

### **PDE7B**

human, recombinant, , N-terminal GST tag  
**Catalog #:** 60071  
**Lot#:** 100127      **Conc.:** 1.5 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0,  
 124 mM NaCl, 2.4 mM KCl, 18 mM  
 glutathione, 3 mM DTT and 10% glycerol.

**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:**

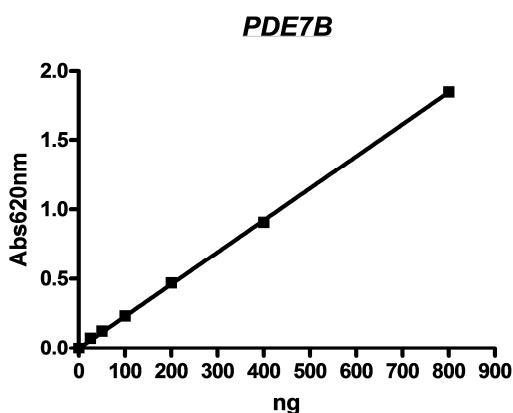
Human PDE7B GenBank Accession No. NM\_018945), (a.a. 107-end) with N-terminal GST tag, MW=66 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

**Specific Activity:** 226 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 Conditions: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl<sub>2</sub>, 0.05% Tween-20, 200 μM cAMP, and PDE7B 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.

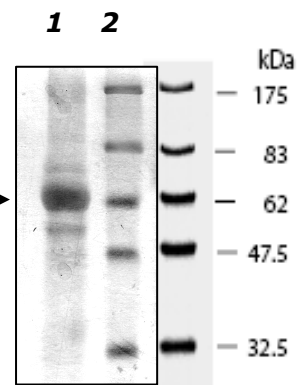
### **Quality Assurance**



**10% SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
 4 μg PDE7B →  
**Lane 2:**  
 Protein Marker  
 BioLabs (#P7708L)

**MW:** 66 kDa.  
**Purity:** >40%



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)