

## Data Sheet

### PDE7A

Rat, recombinant, N-terminal GST-tag

**Catalog #:** 60074

**Lot #:** 101129      **Conc:** 0.34 mg/ml

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

**Stability:** >6 months at -80°C

#### References:

1. Han, P., *et al.* (1998). *Genomics* Mar 1;**48(2)**:275-6.
2. Yang, G., *et al.* (2003). *J. Immunol.* Dec 15;**171(12)**:6414-20.
3. Yamamoto, *et al.* (2006) *Eur. J. Pharmacol.* Jul 10;**541(1-2)**:106-14.

#### Description:

Rat PDE7A (GenBank Accession No. NM\_031080), a.a. 147-482(end), with N-terminal GST-tag, MW=65 kDa, expressed in a Baculovirus-infected Sf9 cell expression system.

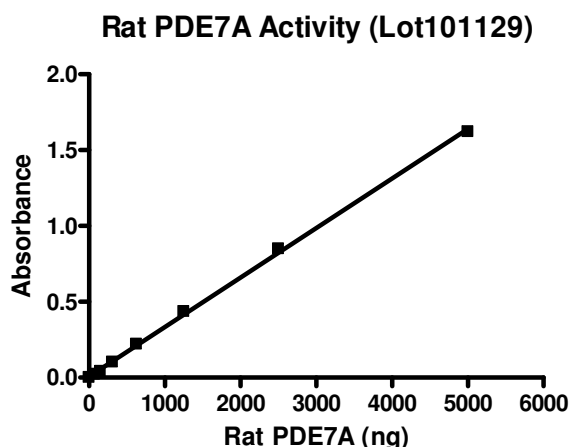
**Specific Activity:** 354 pmol/min/μg.

Assay condition: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl<sub>2</sub>, 0.05% Tween-20, 200 μM cAMP, 0.1 mg/ml BSA, and 0.8-26 ng/μl rat PDE7A 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.

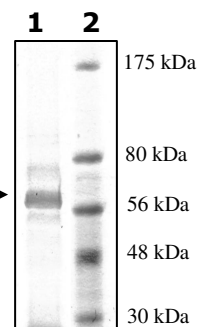
### Quality Assurance



#### 10% SDS-PAGE Coomassie staining

**Lane 1:**  
5μg PDE7A

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



**MW:** 65 kDa.

**Purity:** 30%

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)