



11526 Sorrento Valley Rd SteA2  
San Diego Ca 92121  
Tel: 1.858.202.1401  
1.858.829.3082  
Fax: 1.858.259.5734  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### PDE8A1

human, recombinant, , N-terminal GST tag  
Catalog #: 60080

Formulated in: 25 mM Tris-HCl, pH 8.0,  
100 mM NaCl, 0.05% Tween-20, 50% glycerol,  
and 3 mM DTT.

Stability: >6 months at -80°C

References:

1. Wu P. and Wang P. Proc. Natl. Acad. Sci. U.S.A. 101 (51), 17634-17639 (2004).
2. Gamanuma M. et al., Cell. Signal. 15 (6), 565-574 (2003).

Description:

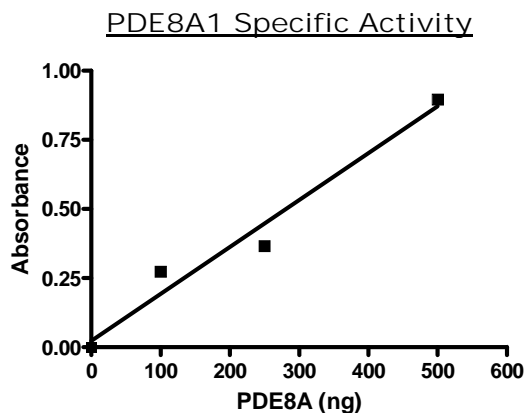
Human PDE8A1 (GenBank Accession No. NM\_002605), full length with N-terminal GST tag, MW= 119 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 437 pmol/min/μg.  
Assay conditions: 10 mM Tris-HCl, pH7.4,  
10 mM MgCl<sub>2</sub>, 1 mM MnCl<sub>2</sub>, 200 μM cAMP, 2.5 kU 5' nucleotidase, 37°C, 20min.

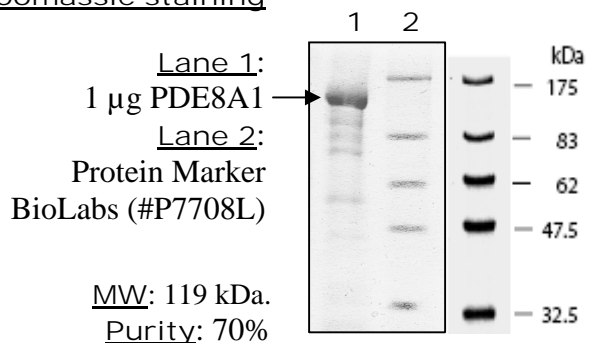
Application:

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

### Quality Assurance



10% SDS-PAGE  
Coomassie staining



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**

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)