



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

Data Sheet

PDE1C

Mouse, recombinant, N-terminal GST tag
Catalog #: 60012

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

Stability: >6 months at -80°C

References:

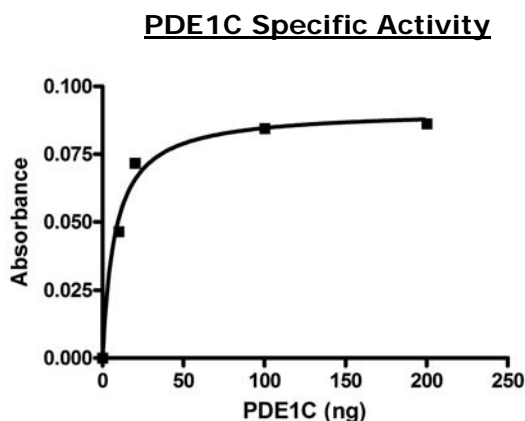
1. Gong S. et al., Nature 425 (6961), 917-925 (2003).
2. Spence S. et al., Biochem. J. 321 (PT 1), 157-163 (1997).

Description: Mouse PDE1C (GenBank Accession No. NM_011054), full length with N-terminal GST tag, MW= 100 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

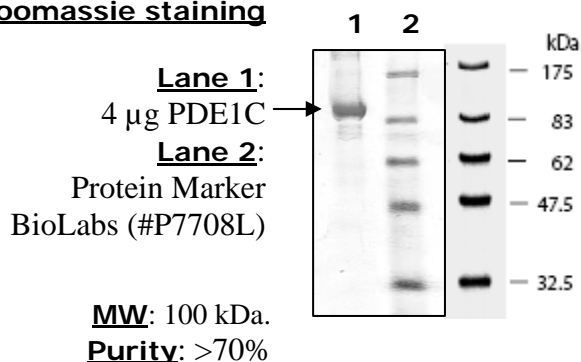
Specific Activity: 5000 U/ug. One unit is defined as the amount of enzyme that will convert 1 pmole of 3', 5'-cAMP to 5' AMP per min at 37 °C in a reaction buffer containing 10 mM Tris-HCl, pH 7.4, 10 mM MgCl₂, 1 mM MnCl₂, 0.4 mM 3', 5'-cAMP and 0.2 ng/ul PDE1C.

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

Please visit our website at: www.bpsbioscience.com