

Data Sheet

PDE6C, GST-Tag (Mouse)

Mouse, Recombinant, N-terminal GST-tag

Catalog #: 60065

Lot #: 161206-2b

Conc.: 0.0063 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCL, 10 mM Glutathione, 3 mM DTT and 20% Glycerol

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary use carrier protein (BSA 0.1 – 0.5%).

References:

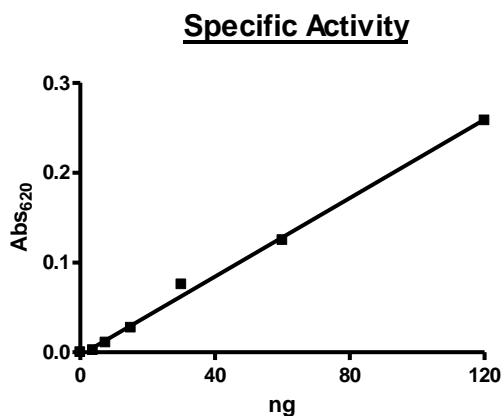
1. Martinez S.E. *et al.*, *J Biol Chem.* 2008 Sep 19; **283(38)**:25913-9.
2. Stearns G. *et al.*, *J Neurosci.* 2007 Dec 12; **27(50)**:13866-74.

Description: Mouse PDE6C, GenBank Accession No. NM_033614, 2-861 (end), with N-terminal GST tag, MW = 126 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: ≥ 600 pmole/min/ μg
 One unit is defined as the amount of enzyme that will convert 1 pmole of 3', 5'-cGMP to 5'-GMP per min at 37°C . Assay condition: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl₂, 0.1 mg/ml BSA, 0.05% Tween-20, 200 μM cGMP, and serial dilutions of PDE6C 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

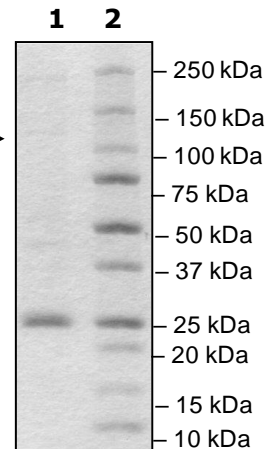


**4-20% SDS-PAGE
 Coomassie staining**

Lane 1: →
 Mouse PDE6C
Lane 2:
 Protein Marker

Purity: $\geq 2\%$

MW: 126 kDa



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
 Please visit our website at: www.bpsbioscience.com