

Data Sheet

PDE11A

Human, recombinant, N-terminal GST tag

Catalog #: 60110

Lot#: 70413

Conc.: 1.7 mg/ml

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

Stability: >6 months at -80°C

References:

1. Loughney, K, *et al.*, *Int. J. Impot. Res.* 2005, **17(4):**320-5.
2. Weeks, JL, *et al.*, *Int. J. Impot. Res.* 2005, **17(1):**5-9.

Description: Human PDE11A (GenBank Accession No. BC112393), full length with N-terminal GST tag, MW= 131 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

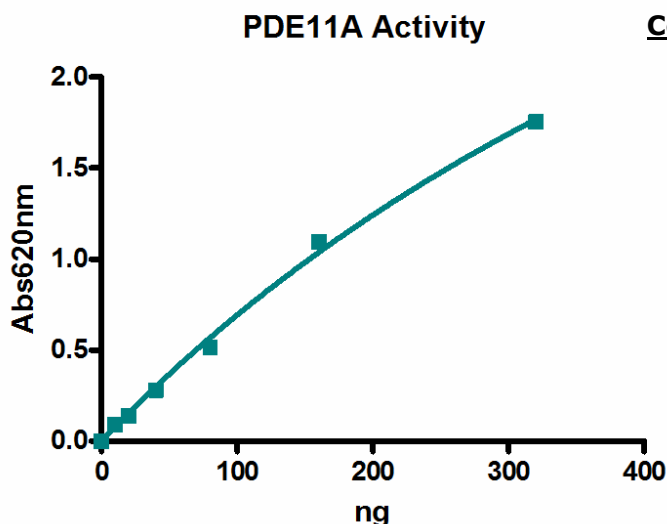
Specific Activity: 1560 pmol/min/μg.

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 condition: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl₂, 0.1 mg/ml BSA, 200 μM cAMP, and PDE11A 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.

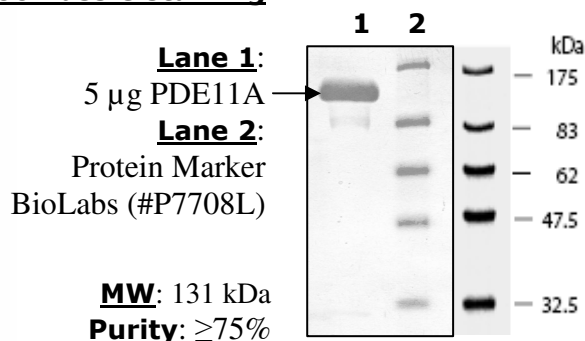
Application:

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

Quality Assurance



**10% SDS-PAGE
 Coomassie staining**



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