

Data Sheet

PDE4D3

Human, recombinant, N-terminal GST tag

Catalog #: 60046

Lot#: 111111GC **Conc.:** 0.42 mg/ml

Formulated in: 50 mM Tris pH 8.0, 138 mM NaCl, 2.7 mM KCl, 0.05% Tween-20, 10% glycerol, and 3mM DTT.

Stability: >6 months at -80°C

References:

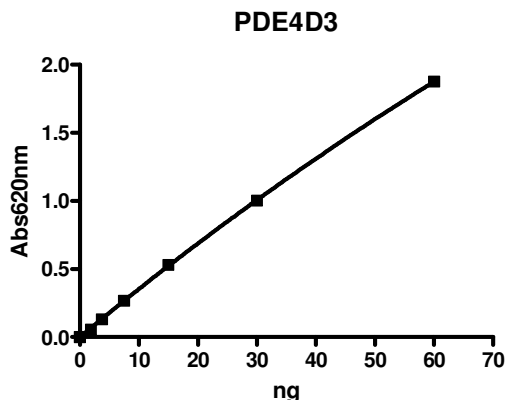
1. Fidani, L., *et al.*, *Eur. J. Neurol.* **14** (7), 745-749 (2007).
2. Baillie, G.S., *et al.*, *Biochem. J.* **404** (1), 71-80 (2007).

Description: Human PDE4D3 (GenBank Accession No. NM_006203), full length, with N-terminal GST tag, MW=102.4 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

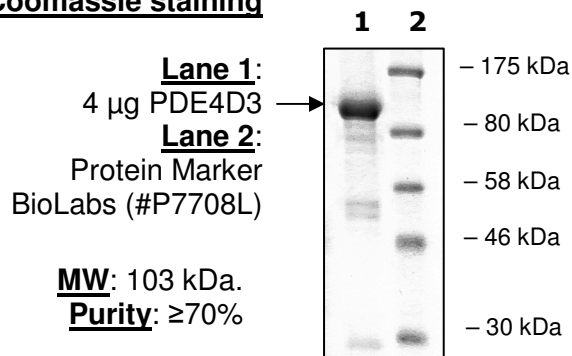
Specific Activity: 5940 pmole/min/μg. Assay condition: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl₂, 1 mM MnCl₂, 200 μM cAMP, 0.1 mg/ml BSA, and 0.2 ng/μl PDE4D3 at 37°C for 20 min. Quantified by 5'-nucleotidase cleaving the 5'-AMP product and releasing the phosphate group which is detected by Malachite Green 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