



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

PDE11A4 (human, recombinant, N-terminal GST tag)
Catalog #: 60110

Description: Human PDE11A4 (GenBank Accession No. BAB62712), full length with N-terminal GST tag, MW= 131 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Source: Baculovirus infected Sf9 cells

Formulated in: 25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20, 50% glycerol, 5 mM glutathione, and 3 mM DTT.

Stability: >6 months at -80°C

Purity: >80% by SDS-PAGE.

Specific Activity: 2008 U/ug. 1 unit is defined as the amount of enzyme that will convert 1 pmole of 3', 5'-cGMP to 5' GMP per min at 30 °C in a reaction buffer of 10 mM Tris-HCl, pH 7.4, 0.2 mM MgCl₂. Assay buffer: 10 mM Tris-HCl, pH 7.4, 1.0 mM MgCl₂, 0.4 mM 3', 5'-cGMP substrate, 1 ng/μl PDE11A4.

Application: useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Reference:

1. Loughney K, Taylor J, and Florio VA *Int J Impot Res.* 2005, 17(4):320-5
2. Weeks JL, Zoraghi R, Beasley A, Sekhar KR, Francis SH, Corbin JD. *Int J Impot Res.* 2005 Jan-Feb;17(1):5-9

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