



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

PDE9A2

Human, recombinant, N-terminal GST tag
Catalog #: 60090

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

Stability: >6 months at -80°C

References:

1. Rentero C. and Puigdomenech P. BMC Mol. Biol. 7, 39 (2006).
2. Huai Q. et al., Journal Proc. Natl. Acad. Sci. U.S.A. 101 (26), 9624-9629 (2004).

Description:

Human PDE9A2 (GenBank Accession No. NM_001001567), full length with N-terminal GST tag, MW=87.7 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 97 pmol/min/μg.
Assay conditions: 10 mM Tris-HCl, pH7.4,
10 mM MgCl₂, 1mM MnCl₂, 400 μM cGMP,
2.5 kU 5' nucleotidase, 5 ng/μl PDE9A,
37°C, 20min.

Application:

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

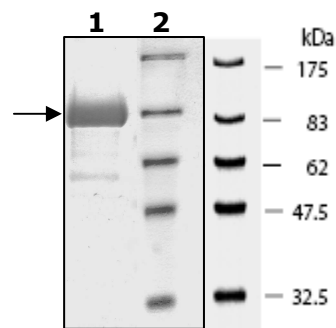
Quality Assurance

**10% SDS-PAGE
Coomassie staining**

Lane 1:
4 μg PDE9A2 →

Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 87.7 kDa.
Purity: >80%



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



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

Assay Protocol

Materials: IMAPTM TR-FRET Screening Express with Progressive Binding Kit from Molecular Devices (R8160); FAM-Cyclic-3',5'-GMP purchased from AXXORA (San Diego, CA); PDE9A2 (BPS Catalog Number 60090).

Methods:

Step 1:

Dilute FAM-cGMP to 200nM in 1X PDE Assay Buffer.
Dilute PDE9A2 enzymes to 0.0125ng/μl in 1X PDE Assay Buffer.

Step 2:

Adding following components to a low binding black plate:
25μl of 200nM FAM-cGMP (Final concentration will be 100nM)
5μl of PDE assay buffer
20μl of PDE9A2 (0.0125ng/μl) (Final amount=0.25ng/reaction)
Mix and incubate at room temperature for 1 hour.

Step 3:

Prepare 1X reagent-binding buffer (75% 1X Binding Buffer A and 25% 1X Binding Buffer B).
Prepare Binding Solution by diluting Binding Reagent with 1X reagent-binding Buffer (1:600).
Add 120μl of Binding Solution to each well and incubate the plate at room temperature for 1 hour.

Step 4:

Measure fluorescence polarization at excitation of 485nm and emission of 520nm in BioTek SynergyTM 2 microplate reader.

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