



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](mailto:info@bpsbioscience.com)

## Data Sheet

### PDE10A1

Human, recombinant, N-terminal GST tag  
Catalog #: 60100

Formulated in: 25 mM Tris-HCl, pH 8.0,  
100 mM NaCl, 0.05% Tween-20, 30% glycerol,  
and 3 mM DTT.

Stability: >6 months at -80°C

References:

1. Gross-Langenhoff, M., et al., J. Biol. Chem. 281 (5), 2841-2846 (2006).
2. Frame, M., et al., Cell. Signal. 13 (10), 735-741 (2001).

Description:

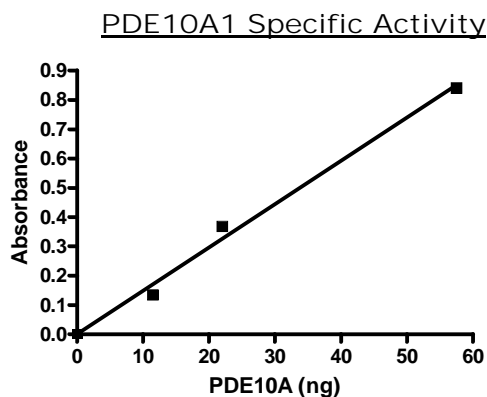
Human PDE10A1 (GenBank Accession No. NM\_006661), full length with N-terminal GST tag, MW= 116 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 2851 pmol/min/μg.  
Assay conditions: 10 mM Tris-HCl, pH7.4,  
10 mM MgCl<sub>2</sub>, 1 mM MnCl<sub>2</sub>, 200 μM cAMP, 2.5 kU 5' nucleotidase, 37°C, 20min.

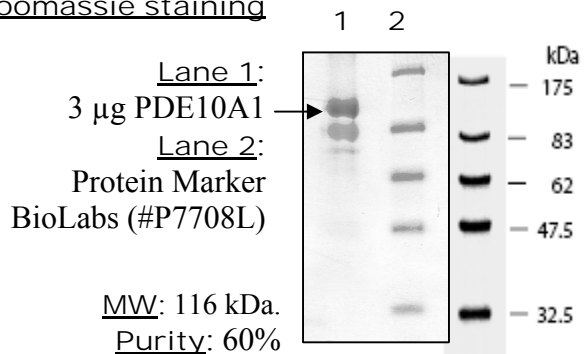
Application:

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 **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://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](mailto:info@bpsbioscience.com)

## Assay Protocol

**Materials:** IMAP<sup>TM</sup> TR-FRET Screening Express with Progressive Binding Kit from Molecular Devices (R8160); FAM-Cyclic-3',5'-AMP from Molecular Devices (R7505); PDE10A1 (BPS Catalog Number 60100).

### **Methods:**

#### **Step 1:**

Dilute FAM-cAMP to 200nM in 1X PDE Assay Buffer.  
Dilute PDE10A1 enzymes to 0.005ng/ $\mu$ l in 1X PDE Assay Buffer.

#### **Step 2:**

Adding following components to a low binding black plate:  
25 $\mu$ l of 200nM FAM-cAMP (Final concentration will be 100nM)  
5 $\mu$ l of 1 x PDE assay buffer  
20 $\mu$ l of PDE10A1 (0.05ng/ $\mu$ l) (Final amount=0.1ng/reaction)  
Mix and incubate at room temperature for 1 hour.

#### **Step 3:**

Prepare 1X reagent-binding buffer (85% 1X Binding Buffer A and 15% 1X Binding Buffer B).  
Prepare Binding Solution by diluting Binding Reagent with 1X reagent-binding Buffer (1:600).  
Add 120 $\mu$ l 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 Synergy<sup>TM</sup> 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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)