

## Data Sheet

### **PDE10A (rat)**

Rat, recombinant, N-terminal GST tag  
 Catalog# 60102  
 Lot#: 80916                      Conc.: 0.81 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
 110 mM NaCl, 2.2 mM KCl, 16 mM  
 glutathione, 20% glycerol, and 3 mM DTT.

**Stability:** >6 months at -80°C

#### **References:**

1. Siuciak, J.A. et al., Neuropharmacology 51 (2), 374-385 (2006).
2. Hu, H. et al., Eur. J. Neurosci. 20 (12), 3351-3363 (2004).

#### **Description:**

Rat PDE10A GenBank Accession No. NM\_022236), amino acids 449-end with N-terminal GST tag, MW=66 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

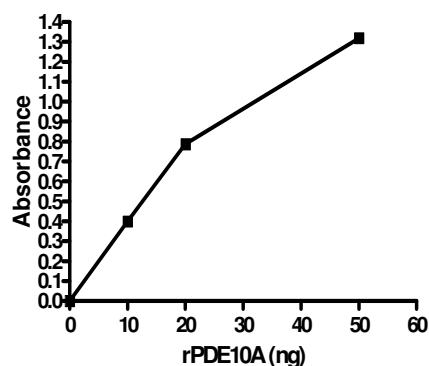
**Specific Activity:** 8339 pmol/min/μg.  
 Assay conditions: 10 mM Tris-HCl, pH7.4, 10 mM MgCl<sub>2</sub>, 1 mM MnCl<sub>2</sub>, 200 μM cAMP, 2.5 kU 5' nucleotidase, 37°C, 20min.

#### **Application:**

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

## **Quality Assurance**

### **PDE10A rat Specific Activity**

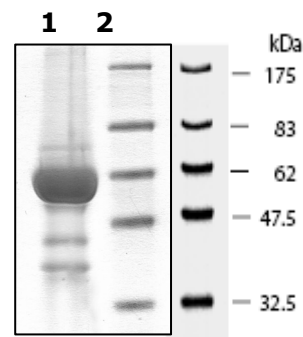


### **10% SDS-PAGE Coomassie staining**

**Lane 1:**  
 7 μg PDE10A rat →

**Lane 2:**  
 Protein Marker  
 BioLabs (#P7708L)

**MW:** 66 kDa.  
**Purity:** >50%



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](mailto:info@bpsbioscience.com)

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