

## Data Sheet

### PDE4A10

Human, recombinant, N-terminal GST tag  
Catalog #: 60038  
Lot#: 120207-GC      Conc.: 0.26 mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 8.0, 300 Glycerol, and 3mM DTT.

**Stability:** >6 months at  $-80^{\circ}\text{C}$

**References:**

1. McCahill, A., et al. *Cell Signal.* 2008. Nov; **20(11)**:2071-83.
2. McPhee, I., et al., *Cell Signal.* 2001. Dec; **13(12)**:911-8.

**Description:** Human PDE4A10, also known as PDE4A protein isoform 6, transcript variant 3 (GenBank Accession No. NM\_001111309), amino acids 2-825(end) with N-terminal GST tag, MW= 117 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

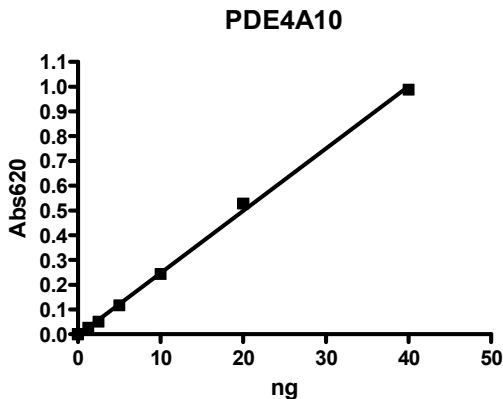
**Specific Activity:** 4362 pmole/min/mg.

One unit is defined as the amount of enzyme that will convert 1 pmole of 3', 5'-cAMP to 5'-AMP per min at  $37^{\circ}\text{C}$ . Assay condition: 10 mM Tris-HCl, pH 7.4, 10 mM  $\text{MgCl}_2$ , 1 mM  $\text{MnCl}_2$ , 200  $\mu\text{M}$  cAMP, and 0.2 ng/ $\mu\text{l}$  PDE4A10 at  $37^{\circ}\text{C}$  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

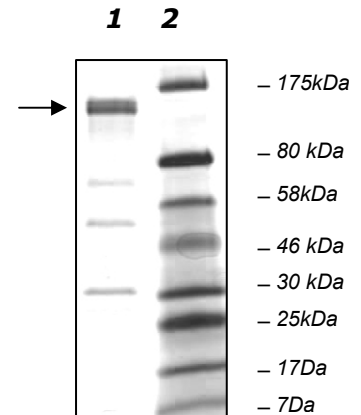


**10% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
1  $\mu\text{g}$  PDE4A10

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

**MW:** 117 kDa  
**Purity:**  $\geq 75\%$



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)