

## PDE12 Protein

Full length recombinant protein expressed in Sf9 cells

**Catalog # P100-30G**

Lot # E117-2

### Product Description

Recombinant full-length human PDE12 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_177966](#).

### Gene Aliases

2'-PDE

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

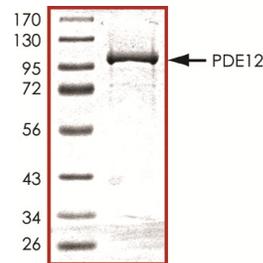
### Scientific Background

PDE12 is 2', 5'-phosphodiesterase 12 protein that belongs to the CCR4/nocturin family. PDE12 cleaves the 2',5'-phosphodiester bond linking adenosines of the 5'-triphosphorylated oligoadenylates. Triphosphorylated oligoadenylates, which are referred to as 2-5A, modulates in the 2-5A system that is one of the major pathways for antiviral and antitumor functions induced by interferons (IFNs) (1). Suppression of PDE12 induces reduction of viral replication in Hela cells, thus counteracting the antiviral pathway probably by inhibiting the 2-5A system.

### References

1. Kubota K. et.al: "Identification of 2'-phosphodiesterase, which plays a role in the 2-5A system regulated by interferon." J. Biol. Chem. 279:37832-37841,2004.

### Purity



The purity was determined to be **>90%** by densitometry. Approx. MW **100kDa**.

## PDE12 Protein

Full length human recombinant protein expressed in Sf9 cells

Catalog Number	P100-30G
Specific Lot Number	E117-2
Purity	>90%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**