

| Catalogue # | Aliquot Size |
|-------------|--------------|
| P89-30H -05 | 5 µg         |
| P89-30H -10 | 10 µg        |
| P89-30H -20 | 20 µg        |

## PDE1A, Active

Full-length recombinant protein expressed in Sf9 cells

**Catalog # P89-30H**

Lot # R148-1

### Product Description

Recombinant full-length human PDE1A was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [NM\\_005019](#).

### Gene Aliases

HCAM1; HSPDE1A; MGC26303

### Formulation

Recombinant protein stored in 50mM MOPS, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . 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.

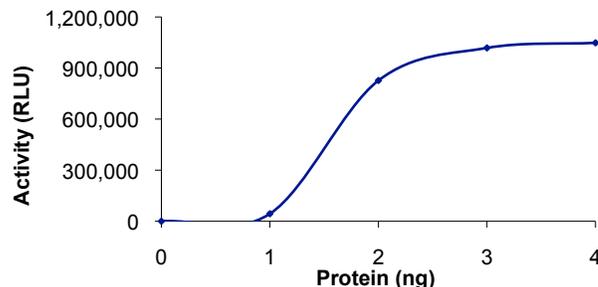
### Scientific Background

PDE1A is a member of the cyclic nucleotide phosphodiesterase family that is stimulated by  $\text{Ca}^{2+}$ -calmodulin. PDE1A is a cGMP-specific phosphodiesterase and hydrolyzes cGMP to 5'-GMP. Calmodulin is tightly bound to PDE1A in human spermatozoa and this results in permanent activation of PDE1A (1). Inhibition of PDE1A function significantly attenuates vascular smooth muscle cell growth by decreasing proliferation via G1 arrest and induces apoptosis (2). In vascular smooth muscle cell, PDE1A is important in regulating growth and survival of these cells and thus can contribute to the neointima formation in atherosclerosis and restenosis.

### References

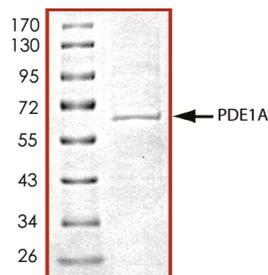
1. Lefièvre L, et al: Presence of cyclic nucleotide phosphodiesterases PDE1A, existing as a stable complex with calmodulin, and PDE3A in human spermatozoa. *Biol Reprod.* 2002 Aug;67(2):423-30.
2. Nagel D J, et al: Role of nuclear  $\text{Ca}^{2+}$ /calmodulin-stimulated phosphodiesterase 1A in vascular smooth muscle cell growth and survival. *Circ Res.* 2006 Mar 31;98(6):777-84.

### Specific Activity



The specific activity of PDE1A was determined to be **12,290 nmol /min/mg** as per activity assay protocol.

### Purity



The purity was determined to be **>90%** by densitometry. Approx. MW **64kDa**.

## PDE1A, Active

Full-length recombinant protein expressed in Sf9 cells

|                     |  |
|---------------------|--|
| Catalog Number      | P89-30H  |
| Specific Activity   | 12,290 nmol/min/mg   |
| Specific Lot Number | R148-1   |
| Purity              | >90%   |
| Concentration       | 0.1µg/µl   |
| Stability           | 1 yr at $-70^{\circ}\text{C}$ from date of shipment  |
| Storage & Shipping  | Store product at $-70^{\circ}\text{C}$ . 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.**

# Activity Assay Protocol

## Reaction Components

### Active PDE1A (Catalog #: P89-30H)

Active PDE1A (0.1µg/µl) diluted with 1X PDE-Glo™ Reaction Buffer and assayed as outlined in sample activity plot. (Note: these are suggested working dilutions and it is recommended that the researcher perform a serial dilution of Active PDE1A for optimal results).

### 100 mM IBMX Solution

Prepare 100 mM of 3-isobutyl-1-methylxanthine (IBMX) in 100% DMSO. Store aliquots at -20°C.

### PDE-Glo™ Phosphodiesterase Assay Kit (Promega, Cat # V1361)

cAMP and cGMP solution, 1 mM  
PDE-Glo™ Reaction Buffer, 5X  
PDE-Glo™ Termination Buffer, 5X  
PDE-Glo™ Detection Buffer, 5X  
Protein Kinase A (PKA)  
Kinase-Glo™ Substrate  
Kinase-Glo™ Buffer

## Assay Protocol

The PDE1A assay is performed using the PDE-Glo™ Phosphodiesterase Assay kit (Promega; Cat# V1361). The assay involves first a PDE1A reaction between an active PDE1A preparation and cyclic nucleotide substrate cGMP. Then PDE-Glo™ Termination Buffer and PDE-Glo™ Detection Buffer (which contains ATP, inactive PKA and PKA substrate) are added to the reaction. The cyclic nucleotide substrate remaining after the PDE1A reaction can bind to the inactive PKA regulatory subunit thereby releasing the active catalytic subunit of PKA. The active catalytic subunit of PKA then catalyzes phosphorylation of the PKA substrate in the presence of ATP which leads to a reduction in ATP level. In the final step, Kinase-Glo™ reagent is added to measure the Luciferase activity towards Luciferin and the luminescent signal produced is related to the amount of ATP remaining which is indirectly related to the activity of PDE1A.

**Step 1.** Thaw the Active PDE1A and PDE-Glo™ Phosphodiesterase Assay Kit reagents on ice.

**Step 2.** Prepare the following working solutions:

- o Diluted active PDE1A with 1X PDE-Glo™ Reaction Buffer on ice
- o 20µM cGMP substrate solution in 1X PDE-Glo™ Reaction Buffer at ambient temperature
- o 1X PDE-Glo™ Termination Buffer in 10 mM IBMX solution at ambient temperature
- o 1X PDE-Glo™ detection solution (mix 8µl PKA with 792µl water and 200µl 5X PDE-Glo™ Detection Buffer). Prepare immediately before use
- o Kinase-Glo™ reagent by adding Kinase-Glo™ Buffer to Kinase-Glo™ Substrate at ambient temperature

**Step 3.** In a polystyrene 96-well plate, add the following components bringing the initial reaction volume up to 25µl:

**Component 1.** 10µl of diluted Active PDE1A (Catalog #P89-30H)

**Component 2.** 2.5µl of 1mM CaCl<sub>2</sub> solution containing 0.15 µg Calmodulin

**Component 3.** 12.5µl of 20µM cGMP solution (0.25 nmol cGMP used per assay)

**Step 4.** Set up a blank control as outlined in step 3 by excluding the addition of the diluted PDE preparation. Replace the PDE preparation with an equal volume of 1X PDE-Glo™ Reaction Buffer.

**Step 5.** Initiate the reaction by adding cGMP substrate solution and incubate the mixture at 30°C for 10 minutes on a plate shaker.

**Step 6.** Terminate the PDE reaction by adding 12.5µl of 1X PDE-Glo™ Termination Buffer. Mix well.

**Step 7.** Add 12.5µl of 1X PDE-Glo™ detection solution. Mix well and then incubate at ambient temperature for 20 minutes.

**Step 8.** After the incubation period, add 50µl of Kinase-Glo™ reagent mix and then incubate at ambient temperature for 10 min.

**Step 9.** Read the polystyrene 96-well reaction plate using the KinaseGlo Luminescence Protocol on a GloMax plate reader (Promega; Cat# E7031).

**Step 10.** Perform a cGMP standard curve. Determine RLU at each concentration. Then calculate the corresponding nmol cGMP remaining after the PDE reaction from the standard curve.

**Step 11.** Calculate the PDE specific activity as outlined below.

### PDE Specific Activity (SA) (nmol/min/mg)

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.**