

PPME1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # P332-30G

Lot # L2110-9

Product Description

Recombinant human mature form of PPME1 (2-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The PPME1 gene accession number is [BC003046](#).

Gene Aliases

PME-1

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM 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.

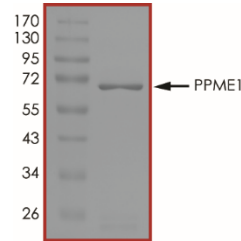
Scientific Background

Protein phosphatase methylesterase 1 (PPME1) conducts the demethylation of the catalytic subunit of protein phosphatase 2A (PP2A) and leads to inactivation of the phosphatase (1). It is demonstrated that PPME1 binds directly to the active site of PP2A and evicts the manganese ions required for the phosphatase activity (2). PPME1 is also reported to protect ERK activity in human malignant glioma (3).

References

- Ogris, E., et al. A protein phosphatase methylesterase (PME-1) is one of several novel proteins stably associating with two inactive mutants of protein phosphatase 2A. *J Biol Chem.* 274(20):14382-91, 1999.
- Xing, Y., et al. Structural mechanism of demethylation and inactivation of protein phosphatase 2A. *Cell.* 133(1):154-63, 2008.
- Puustinen, P., et al. PME-1 protects extracellular signal-regulated kinase pathway activity from protein phosphatase 2A-mediated inactivation in human malignant glioma. *Cancer Res.* 69(7):2870-7, 2009.

Purity



The purity of PPME1 was determined to be **>90%** by densitometry, approx. MW **70 kDa**.

PPME1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	P332-30G
Lot #	L2110-9
Purity	>90%
Concentration	0.1 µg/µl
Stability	1 yr 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
www.signalchem.com

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