

PCMT1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # P314-30G

Lot # L1943-6

Product Description

Recombinant human mature form of PCMT1 (2-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The PCMT1 gene accession number is [BC007501](#).

Gene Aliases

(None)

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.

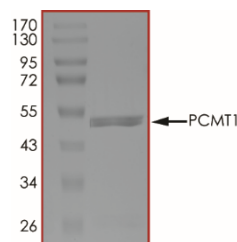
Scientific Background

PCMT1 is a member of the L-isoaspartyl/D-aspartyl protein methyltransferase family, which plays an important role in protein repair (1). Downregulation of PCMT1 leads to hyperactivation of MAPK pathway upon EGF stimulation (2). PCMT1 is regulated by cell-matrix interactions, and linked to various biological processes including wound healing, cell migration, and tumor metastasis (3).

References

1. Tsai, W., et al. Amino acid polymorphisms of the human L-isoaspartyl/D-aspartyl methyltransferase involved in protein repair. *Biochem Biophys Res Commun.* 203(1):491-7, 1994.
2. Kosugi, S., et al. Suppression of protein L-isoaspartyl (d-aspartyl) methyltransferase results in hyperactivation of EGF-stimulated MEK-ERK signaling in cultured mammalian cells. *Biochem Biophys Res Commun.* 371(1):22-7, 2008.
3. Lanthier, J., et al. Regulation of protein L-isoaspartyl methyltransferase by cell-matrix interactions: involvement of integrin alphavbeta3, PI 3-kinase, and the proteasome. *Biochem Cell Biol.* 84(5):684-94, 2006.

Purity



The purity of PCMT1 was determined to be **>95%** by densitometry, approx. MW **51 kDa**.

PCMT1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	P314-30G
Lot #	L1943-6
Purity	>95%
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
www.signalchem.com

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