

PRODUCT INFORMATION



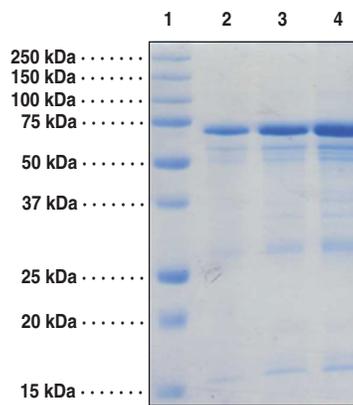
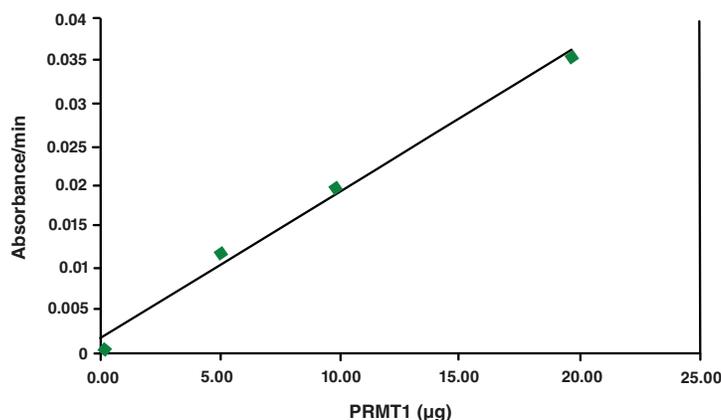
PRMT1 (human recombinant)

Item No. 10350

Overview and Properties

Synonyms: Interferon receptor 1-bound protein 4, Protein Arginine Methyltransferase 1
Source: Active recombinant N-terminal GST-tagged protein expressed in *E. coli*
Amino Acids: 2-371 (full-length)
Uniprot No.: Q99873
Molecular Weight: 69.2 kDa
Storage: -80°C (as supplied); avoid freeze/thaw cycles by aliquoting protein
Stability: ≥6 months
Purity: *batch specific* (≥35% estimated by SDS-PAGE)
Supplied in: *batch specific*
Activity: *batch specific* U/ml one unit is the amount of enzyme required to transfer one methyl group to histone H4 per minute using 100 μmol histone H4 (1-21) peptide at 37°C in Cayman's Methyltransferase Colorimetric Assay Kit (Item No. 700140).
Specific Activity: *batch specific* nanomoles/min/mg Protein
Concentration: *batch specific* mg/ml

Images



Lane 1: MW Markers
Lane 2: PRMT1 (1 μg)
Lane 3: PRMT1 (2 μg)
Lane 4: PRMT1 (4 μg)

Representative gel image shown; actual purity may vary between each batch.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/19/2018

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Protein Arginine Methyltransferases (PRMTs) are S-adenosyl-L-methionine-dependent enzymes that catalyze post-translational modification of arginine residues by methylation in three different ways: monomethylation, symmetric dimethylation, or asymmetric dimethylation at terminal arginine guanidino nitrogen atoms.¹ Class I PRMTs catalyze arginine monomethylation and asymmetric dimethylation. PRMT1 is a class I arginine methyltransferase that methylates arginine residues at a number of glycine and arginine rich regions (GAR motifs) in proteins including histone H4 at arginine 3.²

References

1. Bedford, M.T. Arginine methylation at a glance. *Journal of Cell Science* **120(24)**, 4243-4246 (2007).
2. Wang, H., Huang, Z.Q., Xia, L., *et al.* Methylation of histone H4 at arginine 3 facilitating transcriptional activation by nuclear hormone receptor. *Science* **293**, 853-857 (2001).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM