

## Recombinant PRMT2 protein

**Catalog No:** 31392, 31792

**Expressed In:** Baculovirus

**Quantity:** 20 µg

**Concentration:** 0.4 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant PRMT2 protein is supplied at a concentration of 0.4 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

**Background: PRMT2 (Protein Arginine Methyltransferase 2)** is a member of the protein arginine N-methyltransferase (PRMT) family that is capable of monomethylating and asymmetrically dimethylating arginine residues. PRMT2 methylates the guanidino nitrogens of arginyl residues in proteins such as STAT3, FBL, and histone H4. PRMT2, along with NCOA2, acts as a coactivator of androgen receptor (AR)-mediated transactivation. PRMT2 also acts as a coactivator (with estrogen) of estrogen receptor (ER)-mediated transactivation. PRMT2 enhances PGR, PPAR $\gamma$ , RARA-mediated transactivation and may inhibit NF $\kappa$ B transcription and promote apoptosis. PRMT2 is known to repress E2F1 transcriptional activity in an Rb-dependent manner and may be involved in growth regulation.

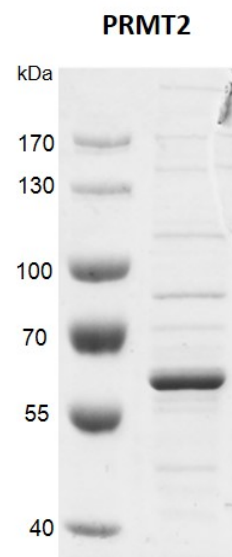
**Protein Details:** Recombinant PRMT2 was expressed in a baculovirus expression system as the full length protein (accession number NP\_996845.1) with an N-terminal FLAG tag. The molecular weight of PRMT2 is 47.6 kDa. The purity of the protein is >75% by SDS-PAGE.

**Application Notes:** Recombinant PRMT2 is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

**Specific Activity:** PRMT2 methylates the guanidino nitrogens of arginyl residues in proteins such as STAT3, FBL, and histone H4. May methylate H3R8 and H4.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



**Recombinant PRMT2 protein gel.** PRMT2 protein was run on an 8% SDS-PAGE gel and stained with Coomassie Blue.