

PRDM2 (KMT8) Protein

Recombinant human protein expressed in Sf9 cells

Catalog # P333-31G

Lot # B1989-5

Product Description

Recombinant PRDM2 (KMT8) (1-340) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The PRDM2 (KMT8) gene accession number is [NM_012231](#).

Gene Aliases

HUMHOXY1; KMT8; MTB-ZF; RIZ; RIZ1; RIZ2

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.

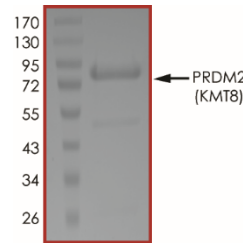
Scientific Background

PRDM2 is a histone methyltransferase that specifically methylates lysine-9 of histone H3. PRDM2 is a retinoblastoma protein-interacting zinc finger protein and a tumor suppressor. Loss of PRDM2 activity through loss of mRNA expression, frameshift, deletion, and missense mutation is commonly found in many types of human cancers (1, 2).

References

1. Du, Y., et al. Hypermethylation in human cancers of the RIZ1 tumor suppressor gene, a member of a histone/protein methyltransferase superfamily. *Cancer Res.* 61(22):8094-9, 2001.
2. Kim, K.C., et al. Inactivation of a histone methyltransferase by mutations in human cancers. *Cancer Res.* 63(22):7619-23, 2003.

Purity



The purity of PRDM2 (KMT8) was determined to be **>85%** by densitometry, approx. MW **80 kDa**.

PRDM2 (KMT8) Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	P333-31G
Lot #	B1989-5
Purity	>85%
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.