
Phospho-PRMT4/CARM1-S228 Polyclonal Antibody

(Catalog #A82997)

Background

This gene belongs to the protein arginine methyltransferase (PRMT) family. The encoded enzyme catalyzes the methylation of guanidino nitrogens of arginyl residues of proteins. The enzyme acts specifically on histones and other chromatin-associated proteins and is involved in regulation of gene expression. The enzyme may act in association with other proteins or within multi-protein complexes and may play a role in cell type-specific functions and cell lineage specification. A related pseudogene is located on chromosome 9.

Description

Phospho-PRMT4/CARM1-S228 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thimerosal, 50% glycerol, pH7.3.

Specificity

Human

Isotype

IgG

Uniprot ID

Q86X55

Purification

Affinity Purified

Immunogen

Synthetic peptide. Please contact EpigenTek for more information.

Storage

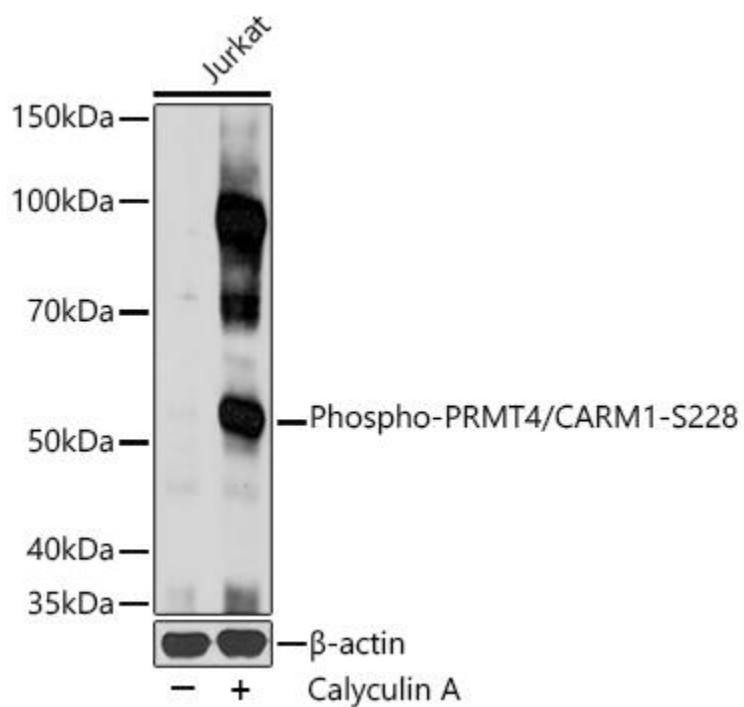
Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

PRMT4; Phospho-PRMT4/CARM1-S228

Application

WB, ELISA; Recommended dilution: WB, 1:500 - 1:1000



Western blot analysis of lysates from Jurkat cells, using Phospho-PRMT4/CARM1-S228 Polyclonal Antibody at 1:1000 dilution. Jurkat cells were treated with Calyculin A (100 nM) at 37°C for 30 minutes.

Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25µg per lane.

Blocking buffer: 3% nonfat dry milk in TBST.

Exposure time: 10s.