

## CARM1 antibody (pAb)

**Catalog Nos:** 39251, 39252

**RRID:** AB\_2793205

**Isotype:** Serum

**Application(s):** ChIP, IHC, IP, WB

**Reactivity:** Human, Mouse, Rat

**Volumes:** 200 µl, 10 µl

**Purification:** None

**Host:** Rabbit

**Molecular Weight:** 60 kDa

**Background:** CARM1 (Coactivator-associated arginine methyltransferase 1 or CARM 1) is a protein arginine N-methyltransferase that catalyzes the mono- and asymmetric dimethylation of arginine residues in its substrate proteins. It serves as a coactivator of transcription and also plays a role in nuclear receptors. CARM1 methylates histone H3 at Arg2, Arg17 and Arg26; it also methylates a number of non-histone substrates involved in the regulation and mechanism of gene expression.

**Immunogen:** This CARM1 antibody was raised against a peptide corresponding to the C-terminus of mouse CARM1.

**Buffer:** Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

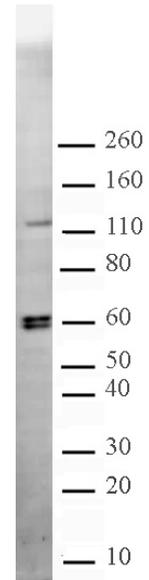
Applications Validated by Active Motif:

WB: 1:250 - 1:1,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

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



**CARM1 pAb tested by Western blot.**  
Whole-cell extract of HEK293 cells (10 µg) probed with CARM1 pAb at a dilution of 1:500.