

## PRDM1 / BLIMP1 antibody (mAb)

---

**Catalog Nos:** 61167, 61168

**RRID:** AB\_2615011

**Clone:** ROS195

**Isotype:** IgG1

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

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 77 & 95 kDa

---

**Background:** PRDM1 / BLIMP1 is a transcriptional repressor that binds to specific target sites in DNA and recruits co-repressors. PRDM1 / BLIMP1 is a master regulator for B cell terminal differentiation, is essential for normal development and immunity and a putative tumor suppressor in diffuse large B cell lymphomas (DLBCLs). PRDM1 / BLIMP1 specifies primordial germ cells through the silencing of the default somatic pathway. PRDM1 / BLIMP1 also acts as a master regulator of intestinal maturation.

**Immunogen:** This PRDM1 / BLIMP1 antibody was raised against full-length recombinant human PRDM1 / BLIMP1 protein.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

**Application Notes:**

**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.