

## PML antibody (mAb)

---

**Catalog Nos:** 39773, 39774

**RRID:** AB\_2793337

**Clone:** 36.1-104

**Isotype:** IgG

**Application(s):** ICC, IF, IP, WB

**Reactivity:** Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 110 kDa

---

**Background:** PML (Promyelocytic Leukemia) protein is a transcription factor originally identified due to a chromosomal aberration fusing it to the Retinoic Acid Receptor alpha (RAR-A) in cases of Acute Promyelocytic Leukemia (APL). Besides transcriptional regulation, PML is involved in many processes, including the response to DNA damage, cell cycle regulation, viral infection response and apoptosis. These functions of PML involve the localization of PML to unique sub-nuclear structures called PML Nuclear Bodies (PML NBs). Over fifty proteins are known to be localized to PML NBs, either transiently or constitutively. These Nuclear Bodies are disrupted in PML knockout mice. PML is subject to a number of post-translational modifications, including phosphorylation, ubiquitylation and sumoylation.

**Immunogen:** This PML antibody was raised against a recombinant protein corresponding to amino acids 1-581 of mouse PML.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 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.