

## ING1 antibody (mAb)

---

**Catalog Nos:** 61459, 61460

**RRID:** AB\_2793646

**Clone:** Cab1

**Isotype:** IgG2a

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

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein A Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 47 kDa

---

**Background:** **ING1** (Inhibitor Of Growth Family, Member 1) is a member of the inhibitor of growth (ING) family of proteins that play key roles in cell cycle arrest, apoptosis, cell aging, and the DNA damage response. ING1 contains a PHD domain that interacts with Histone H3K4me3. ING1 cooperates with p53/TP53 in the negative regulatory pathway of cell growth by modulating p53-dependent transcriptional activation. Implicated as a tumor suppressor.

**Immunogen:** This antibody was raised against a recombinant protein corresponding to amino acids 67 - 279 of human ING1.

**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.