

## Ago3 antibody (mAb)

---

**Catalog Nos:** 39787, 39788

**RRID:** AB\_2793344

**Clone:** 4B1-F6

**Isotype:** IgG2a

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

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 110 kDa

---

**Background:** Argonaute 3 (Ago3, aka EIF2C3) is a member of the Argonaute family of proteins involved in RNAi mediated gene silencing through siRNA and miRNA effectors. The Argonaute proteins are part of RISC, the RNAi Induced Silencing Complex. Ago3, however, lacks RNase activity and does not appear to cleave target mRNA molecules. Ago3 has both a PAZ and a PIWI domain, indicative of the Argonaute family of proteins.

**Immunogen:** This Ago3 (Argonaute 3) antibody was raised against a recombinant protein corresponding to amino acids 1-175 of human Ago3.

**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.