

Ago1 antibody (mAb)

Catalog Nos: 61071, 61072

RRID: AB_2793495

Clone: 4B8

Isotype: IgG2a

Application(s): IP, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 100 kDa

Background: The **Argonaute** family proteins (**Ago1**, **Ago2**, **Ago3** and **Ago4**) are involved in **RNAi** mediated gene silencing through siRNA and miRNA effectors. The Argonaute proteins are part of **RISC**, the RNAi Induced Silencing Complex. Binding of RISC can inhibit mRNA translation and results in the degradation of the mRNA through the action of **Ago2**-mediated target cleavage. **Ago1** is non-essential for siRNA-directed target RNA cleavage but is needed for the production of mature miRNA that affects miRNA-directed RNA cleavage. Argonaute family proteins contain a **PAZ** and a PIWI domain.

Immunogen: This Ago1 antibody was raised against a peptide from the N-terminus of human Argonaute 1 (Ago1) protein.

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.