

## Ago1/2/3 antibody (mAb)

**Catalog Nos:** 39937, 39938

**RRID:** AB\_2793399

**Clone:** 2A8

**Isotype:** IgG

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

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 95 kDa

**Background:** The Argonaute family of 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. Ago3 lacks RNase activity and does not appear to cleave target mRNA molecules. Argonaute family proteins contain a PAZ and a PIWI domain.

**Immunogen:** This Ago1/2/3 antibody was raised against a recombinant protein corresponding to amino acids 47-859 of human Ago2.

**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:

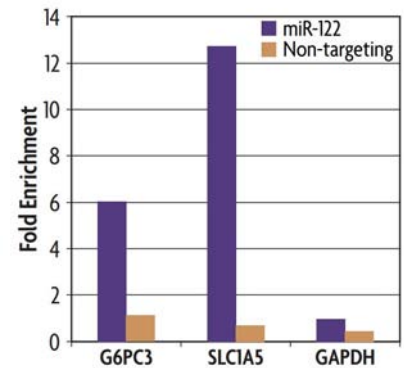
Applications Validated by Active Motif:

IP: 5 µg per IP

This antibody is also available as an AbFlex<sup>®</sup> engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91217.

**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.



**Ago1/2/3 antibody (mAb) (Clone 2A8) tested by immunoprecipitation.** The Active Motif miRNA IP Kit was used on samples of HT1080 cells that had been transfected with a miR-122 or a non-targeting miRNA control for 8 hours. Following IP using the Ago1/2/3 antibody or negative control IgG, qRT-PCR was performed on the samples using primers for G6PC3 and SLC1A5, which are known targets of miR-122, and for GAPDH, a common housekeeping gene that is not known to be targeted by miR-122.