

Ago2 antibody (mAb)

Catalog Nos: 39853, 39854

RRID: AB_2793369

Clone: 1B1-E2H5

Isotype: IgG2a

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

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 100 kDa

Background: Argonaute 2 (Ago2, aka EIF2C2) 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. Binding of RISC can inhibit mRNA translation and results in the degradation of the mRNA through the action of Ago2-mediated target cleavage.

Immunogen: This Ago2 antibody was raised against a recombinant protein corresponding to amino acids 1-148 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:

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.