

Dicer antibody (mAb)

Catalog Nos: 39817, 39818

RRID: AB_2793353

Clone: N167/7

Isotype: IgG1

Application(s): WB

Reactivity: Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 245 kDa

Background: Dicer is a double-stranded RNase and part of the micro-RNA (miRNA) loading complex (miRLC), which is composed of Dicer, Ago2 and TARBP2. Dicer is required for formation of the RNA induced silencing complex (RISC). Dicer and TARBP2 are required to process precursor miRNAs (pre-miRNAs) to mature miRNAs and then load them onto Ago2. Dicer also cleaves double-stranded RNA to produce short interfering RNAs (siRNAs) which target the destruction of complementary RNAs.

Immunogen: This Dicer antibody was raised against a recombinant protein corresponding to amino acids 1638-1899 of mouse Dicer.

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:

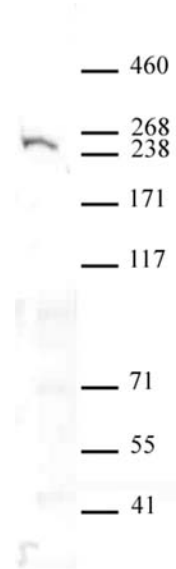
Applications Validated by Active Motif:

WB: 1 - 2 µg/ml dilution

This antibody is also available as an AbFlex® engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91241.

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.



Dicer antibody (mAb) tested by Western blot.
Cytoplasmic extract of mouse ES cells (40 µg per lane) probed with Dicer antibody (2 µg/ml).