

AbFlex[®] Dicer antibody (rAb)

Catalog Nos: 91241, 91242

RRID: AB_2793815

Isotype: IgG2a

Application(s): WB

Reactivity: Mouse

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 240 kDa

Background: AbFlex[®] antibodies are recombinant antibodies (rAbs) that have been generated using defined DNA sequences to produce highly specific, reproducible antibodies. Each AbFlex antibody contains a 6xHis Tag, a Biotinylation Tag for enzymatic biotin conjugation using the biotin ligase, BirA, and a sortase recognition motif (LPXTG) to attach a variety of labels directly to the antibody including fluorophores, enzymatic substrates (HRP, AP), peptides, drugs as well as solid supports. AbFlex[®] Dicer antibody was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells.

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 antibody was raised against amino acids 1638-1899 of mouse Dicer (aka dicer 1, ribonuclease type III).

Buffer: Purified IgG in 140 mM Hepes, pH 7.5, 70 mM NaCl, 32 mM NaOAc, 0.035% sodium azide, 30% glycerol. Sodium azide is highly toxic.

Application Notes:

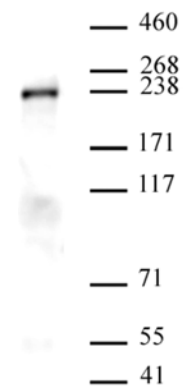
Applications Validated by Active Motif:

WB: 0.5 - 2 µg/ml

AbFlex[®] recombinant antibodies are genetically derived from DNA sequences of parental hybridoma clones. For details on the parental clone, see Catalog No. 39817.

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.



AbFlex[®] Dicer antibody (rAb) tested by Western blot. 40 µg of mouse embryonic stem cell extract was run on SDS-PAGE and probed with AbFlex Dicer antibody at 2 µg/ml.