

AbFlex[®] Mili / PiwiL2 antibody (rAb)

Catalog Nos: 91253, 91254

RRID: AB_2793820

Isotype: IgG2a

Application(s): WB

Reactivity: Mouse

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 110 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[®] Mili / PiwiL2 antibody was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells.

Mili / PiwiL2 is a member of the PIWI/AGO subfamily of RNA-binding proteins. **Mili / PiwiL2** contains a Piwi and an AGO domain and binds a class of RNA called piwi-interacting RNA (piRNA), which are 24 to 30 nucleotides in length and are derived primarily from transposons and other repeated sequence elements. **Mili / PiwiL2** is required for methylation and inactivation of retrotransposons. Mice deficient in Mili / PiwiL2 are sterile, as it is also required for spermatogenesis. It also positively regulates translation, and is required for germ line stem cell division and differentiation.

Immunogen: This antibody was raised against full length mouse MILI / PIWIL2 (Mili like / piwi-like 2) recombinant protein.

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:

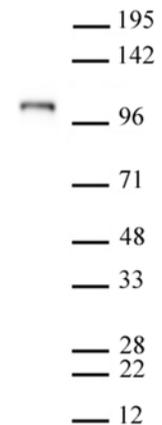
Applications Validated by Active Motif:

WB: 0.2 - 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. 61143.

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[®] Mili / PiwiL2 antibody (rAb) tested by Western blot.

20 µg of mouse testis tissue extract was run on SDS-PAGE and probed with AbFlex Mili / PiwiL2 antibody at 2 µg/ml.