

## Mili / PiwiL2 antibody (mAb)

---

**Catalog Nos:** 61143, 61144

**RRID:** AB\_2793524

**Clone:** 17.8

**Isotype:** IgG1

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

**Reactivity:** Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 97 kDa

---

**Background:** 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 derived 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 Mili / PiwiL2 antibody was raised against full-length recombinant mouse MILI / PIWIL2 protein.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

**Application Notes:**

This antibody is also available as an AbFlex® engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91253.

**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.