

## EED antibody (mAb)

**Catalog Nos:** 61203, 61204

**RRID:** AB\_2615071

**Clone:** 41D

**Isotype:** IgG2a

**Application(s):** ChIP, ChIP-Seq, IHC, WB

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein A Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 57 kDa

**Background:** EED (extraembryonic ectoderm) is a component of the Polycomb group (PcG) multiprotein complex PRC2. The PRC2 complex also contains EZH2 and SUZ12 and is recruited to Polycomb response elements and results in the methylation of histone H3 and lysine 27 and the silencing of gene expression. PRC2 initiates repression of the Hox genes during development and is involved in the maintenance of stem cell pluripotency. EED is a member of the superfamily of WD-40 repeat protein family and is required for PRC2-mediated H3K27 methylation. The PRC2 complex may also recruit platform DNMTs, thereby linking two systems of epigenetic repression.

**Immunogen:** This EED antibody was raised against full-length recombinant human EED protein.

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

### Application Notes:

Applications Validated by Active Motif:

ChIP: 4 µl per ChIP

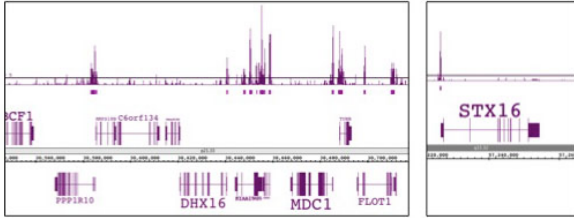
ChIP-Seq: 4 µl each

IHC: 2 - 8 µg/ml dilution

This antibody is also available as an AbFlex<sup>®</sup> engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91135.

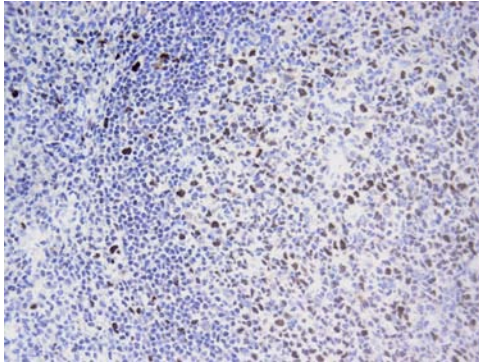
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This 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.



#### EED antibody (mAb) tested by ChIP-Seq.

ChIP was performed using the ChIP-IT® High Sensitivity Kit (Cat. No. 53040) with chromatin from a human B cell lymphoma cell line (4.5 million cells) and 4 µl of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 32 million sequence tags were mapped to identify EED binding sites. The image on the left shows EED binding across a 1.6 million bp region on chromosome 6. The image on the right shows EED binding at the STX16 start site.



#### EED mAb tested by Immunohistochemistry .

EED detection by Immunohistochemistry. The analysis was performed using human tonsil paraffin section and the EED mAb.