

## AbFlex® EED antibody (rAb)

**Catalog Nos:** 91135, 91136

**RRID:** AB\_2793786

**Application(s):** ChIP, ChIP-Seq, ELISA

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Ni-NTA

**Host:** Mouse

**Isotype:** IgG2a

**Concentration:** 1 µg/µl

**Molecular Weight:** 57 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® EED antibody was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells.

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 140 mM Hepes, pH 7.5, 70 mM NaCl, 32 mM NaOAc, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

Applications Validated by Active Motif:

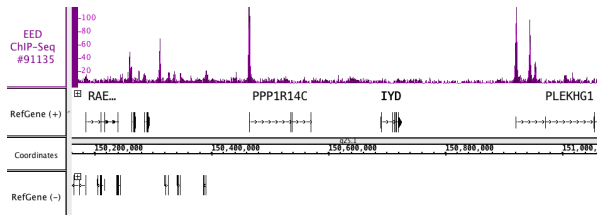
ChIP-Seq: 4 µg per ChIP

Bead-based ELISA: 4 - 26 ng/ml

AbFlex® recombinant antibodies are genetically derived from DNA sequences of parental hybridoma clones. For details on the parental clone, see Catalog No. 61203.

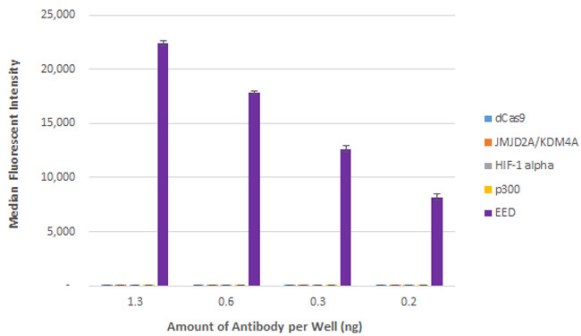
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This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

ChIP was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with 30  $\mu$ g chromatin from an acute large B-cell lymphoma cell line and 4  $\mu$ g of EED antibody. ChIP DNA was sequenced on the Illumina HiSeq and 25 million sequence tags were mapped to identify EED binding sites. The image shows binding across a region of chromosome 6.



AbFlex<sup>®</sup> EED antibody (rAb) tested by Luminex bead-based specificity analysis.

Luminex bead-based specificity analysis was used to confirm the specificity of AbFlex<sup>TM</sup> EED antibody (rAb) for EED. Various proteins were conjugated to MagPlex Luminex beads and incubated with various amounts of AbFlex<sup>TM</sup> EED antibody (rAb). Protein-bound antibody was detected with anti-mouse IgG-Phycoerythrin and read in a Luminex instrument.

Luminex<sup>®</sup> is a registered trademark of Luminex Corporation.