

TAL-1 antibody (mAb)

Catalog Nos: 61259, 61260

RRID: AB_2793572

Clone: BTL73

Isotype: IgG1

Application(s): EMSA, IHC, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 42 kDa

Background: The TAL-1 (T-cell acute lymphocytic leukemia 1, BHLHA17, SCL, TCL5) gene encodes a protein that is implicated in the genesis of hematopoietic malignancies. It may play an important role in hematopoietic differentiation and serves as a positive regulator of erythroid differentiation. TAL-1 is a member of the helix-loop-helix family of transcription factors. Overexpression of TAL-1 is the most common molecular abnormality found in human T cell leukemia.

Immunogen: This TAL-1 antibody was raised against full-length recombinant human TAL-1 protein.

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

Application Notes:

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.