

Recombinant MLLT1 / ENL (1-148) protein

Catalog No: 81098, 81898

Lot No: 34217001

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 0.75 µg/µl

Source: Human

Buffer Contents: Recombinant MLLT1 / ENL (1-148) protein is supplied in 25 mM Tris- HCl pH 8.0, 300 mM NaCl, 10% glycerol and 0.5 mM TCEP.

Background: MLLT1 / ENL (Myeloid/Lymphoid Or Mixed-Lineage Leukemia; Translocated To, 1), also known as YEATS1 or LTG19, is a component of the super elongation complex (SEC), a complex required to increase the catalytic rate of RNA polymerase II transcription by suppressing transient pausing by the polymerase at multiple sites along the DNA.

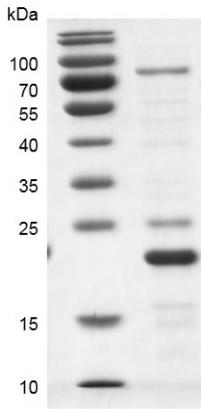
Protein Details: Recombinant human MLLT1 / ENL (1-148) that includes amino acids 1-148 (accession number NP_005925.2) was expressed in *E. coli* cells with an N-terminal 6×His tag and a C-terminal Flag tag. The molecular weight of the protein is 21.9 kDa.

Application Notes: Recombinant MLLT1 / ENL (1-148) protein is suitable for use in enzyme kinetics, inhibitor screening, and selectivity profiling.

Binding Assay Conditions: 3 µM H3K27cr (crotonylated Lysine 27) peptide was incubated with different concentrations of MLLT1 / ENL (1-148) protein in 10 µl reaction system containing 50mM HEPES-NaOH pH7.5, 0.1% BSA for 1 hour, then 10 µl FLAG antibody and SA-XL665 mixture (each 1:100 dilution in Binding Buffer) was added to each reaction system and incubated for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

MLLT1 / ENL (1-148)



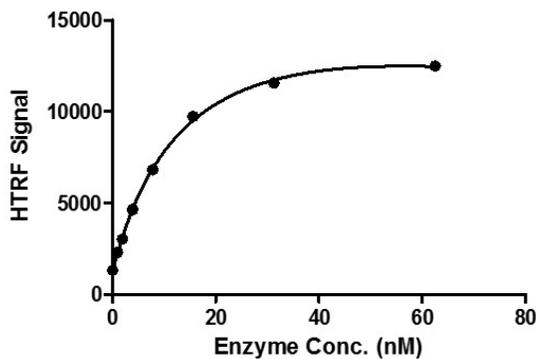
Recombinant MLLT1 / ENL (1-148) protein gel

12.5% SDS-PAGE Coomassie staining

MW: 21.9 kDa

Purity: >70%

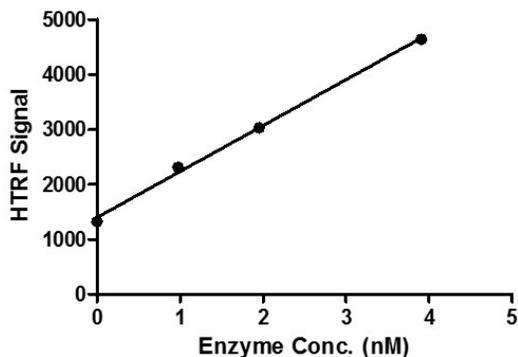
MLLT1 / ENL (1-148) Titration



Recombinant MLLT1 / ENL (1-148) activity assay

H3K27cr (crotonylated Lysine 27) peptide was incubated with different concentrations of MLLT1 / ENL (1-148) protein in reaction buffer for 1 hour, then FLAG antibody and SA-XL665 mixture was added to each reaction system for 30 min. All operations and reactions were performed at room temperature. HTRF assay was used for detection.

MLLT1 / ENL (1-148) Titration



Recombinant MLLT1 / ENL (1-148) activity assay

H3K27cr (crotonylated Lysine 27) peptide was incubated with different concentrations of MLLT1 / ENL (1-148) protein in reaction buffer for 1 hour, then FLAG antibody and SA-XL665 mixture was added to each reaction system for 30 min. All operations and reactions were performed at room temperature. HTRF assay was used for detection.