

ENL YEATS Domain, Y78A, Recombinant Human



EpiCypher®

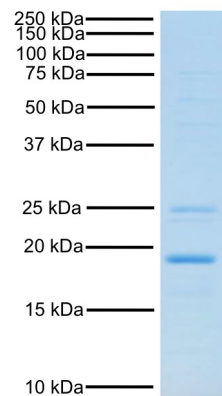
Catalog No 15-0070
Lot No 172562001
Pack Size 100 µg

Type YEATS **Expressed In** *E. coli*

Mol. Wgt. 19 kDa **Epitope Tag** 6xHis

Product Description:

ENL YEATS Domain, Y78A, Recombinant Human (MLLT1, LTG19, YEATS1, accession Q03111, amino acids 1 to 150) containing an N-terminal 6xHis tag, expressed in *E. coli*. The ENL protein is part of the super elongation complex regulating RNA Pol II transcription. A chromosomal translocation resulting in an MLL-ENL fusion is involved in pediatric Acute Lymphoblastic Leukemia. The Y78A substitution ablates ENL YEATS domain binding activity.



Protein Gel Data: ENL YEATS Domain, Y78A, Recombinant Human (1 µg) was run on an SDS- PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.

Formulation:

Recombinant 6xHis-tagged protein at 1 mg/mL in 50 mM Tris-HCl pH 7.5, 200 mM NaCl, 20% glycerol

Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.

Application Notes:

ENL YEATS Domain, Y78A, Recombinant Human is a useful negative control for protein binding and screening experiments with ENL YEATS Domain (EpiCypher 15-0069) with acetylated

References:

This product is for *in vitro* research use only and is not intended for use in humans or animals.