

# ENL YEATS Domain, Recombinant Human



## EpiCypher®

**Catalog No** 15-0069  
**Lot No** 21209006-03  
**Pack Size** 100 µg

**Type** YEATS  
**Expressed In** *E. coli*  
**Mol. Wgt.** 19 kDa  
**Epitope Tag** 6xHis

### Product Description:

ENL YEATS Domain (MLLT1, LTG19, YEATS1, accession Q03111, amino acids 1 to 150) containing a C-terminal 6xHis tag, expressed in *E. coli*. The ENL protein is part of a super elongation complex (SEC) regulating RNA Pol II transcription, and a chromosomal translocation resulting in an MLL-ENL fusion is involved in pediatric Acute Lymphoblastic Leukemia [1]. ENL YEATS Domain preferentially interacts with histone H3 crotonylated lysines over acetylated lysines [2].

### Formulation:

Recombinant 6xHis-tagged protein at 1 mg/mL in 20 mM Tris HCl pH 7.5, 500 mM NaCl, 5% glycerol, 1 mM DTT. Molarity = 52.6 µM

### Storage and Stability:

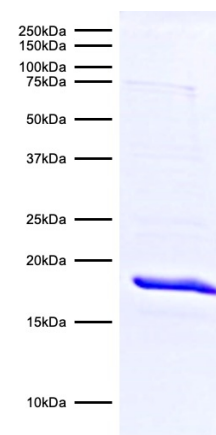
Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

### Application Notes:

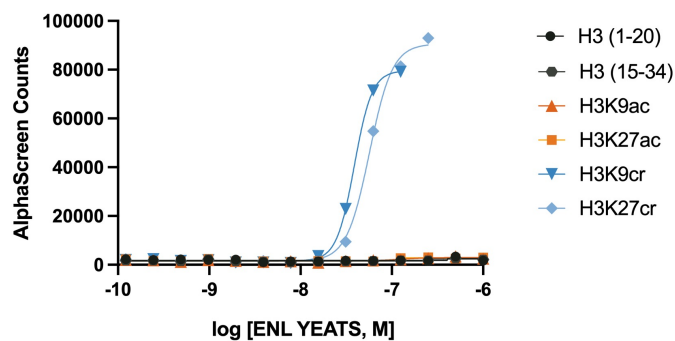
ENL YEATS Domain is useful for protein binding and screening experiments examining crotonylated and acetylated protein substrates.

### References:

- [1] Lin C et al. (2010) *Mol Cell* 37: 429-437.  
[2] Sabari et al. (2016) *Nat Rev Mol Cell Biol* 18: 90-101.



**Protein Gel Data:** ENL YEATS Domain (1 µg) was resolved via SDS-PAGE gel and stained with Coomassie blue to demonstrate the purity of the protein. The migration and molecular weight of the protein standards are indicated.



**Protein Interaction Data:** ENL YEATS Domain shows preferential binding to crotonylated H3 peptides when assayed using AlphaScreen®.

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