

AF9 YEATS Domain, Recombinant Human



EpiCypher®

Catalog No 15-0071
Lot No 21190002-01
Pack Size 100 µg

Type YEATS
Expressed In *E. coli*
Mol. Wgt. 19.6 kDa
Epitope Tag His

Product Description:

AF9 YEATS Domain (MLLT3, YEATS3, accession P42568, amino acids 1 to 149) expressed in *E. coli* containing an N-terminal 6xHis tag. The AF9 protein is part of the super elongation complex and also associates with the histone methyltransferase DOT1L [1]. Translocation of the AF9 gene to the MLL gene, resulting in an MLL/AF9 fusion, is the most common chromosomal rearrangement involving MLL in de novo AML [2].

Formulation:

Recombinant 6xHis-tagged protein at 0.928 mg/mL in 20mM Tris HCl pH 7.5, 500 mM NaCl, 5% glycerol, 1 mM DTT. Molarity = 47.35 µM.

Storage and Stability:

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

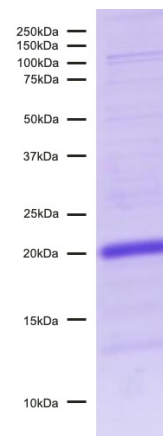
Application Notes:

AF9 YEATS Domain is useful for protein binding and screening experiments examining crotonylated and acetylated protein substrates.

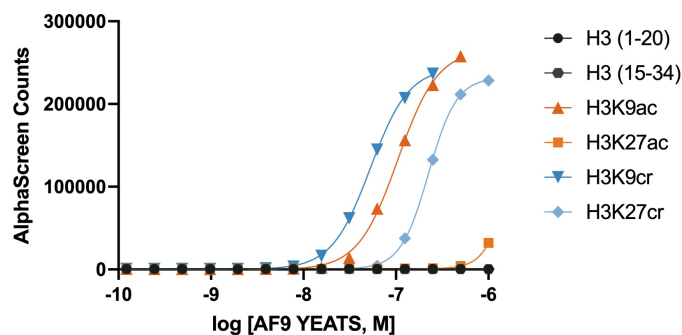
References:

- [1] Li Y et al. (2014) *Cell* 159: 558-571.
- [2] Krivtsov AV et al. (2006) *Nature* 422: 818-822

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



Protein Gel Data: AF9 YEATS Domain (1 µg) was resolved via PAGE gel and stained with Coomassie blue to demonstrate the size and purity of the protein. The migration and molecular weight of the protein standards are indicated.



Protein Interaction Data: AF9 YEATS Domain demonstrates preferential binding to H3K9cr, H3K9ac, and to a lesser extent H3K27cr peptides, but no significant binding to H3K27ac when analyzed by AlphaScreen®.