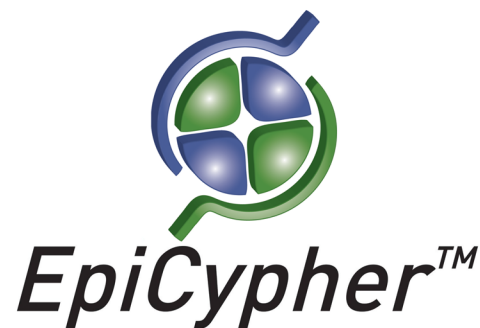


AF9 YEATS Domain, Recombinant Human

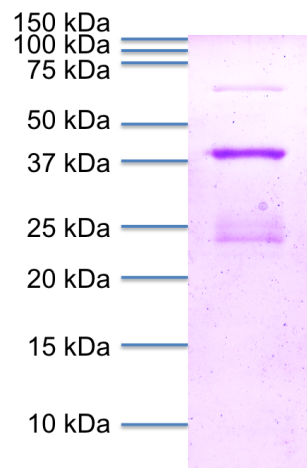
Catalog No. 15-0045
Lot No. 14336001
Pack Size 100 µg



Type YEATS **Expressed In** *E. coli*
Mol. Wgt. 41 kDa **Epitope Tag** GST

Product Description:

Recombinant human AF9 (MLLT3, YEATS3, accession P42568, amino acids 1 to 149) containing an N-terminal GST tag, expressed in *E. coli*. The AF9 protein is part of the super elongation complex and also associates with the histone methyltransferase DOT1L. 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. The YEATS domain of binds strongly to histone H3K9 acetylation and, to a lesser extent, H3K27 and H3K18 acetylation.



Protein Gel Data: AF9 YEATS Domain, Recombinant Human (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.

Formulation:

Recombinant protein at 1 mg/ml in 20 mM Tris-HCl pH 7.5, 150 mM NaCl, 1mM EDTA, 1 mM DTT, 20% glycerol.

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, Recombinant Human is useful for protein binding and screening experiments examining acetylated protein substrates.

References:

Li Y et al (2013) *Cell* 159: 558-571.

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