

YEATS2 YEATS Domain, Recombinant Human



EpiCypher®

Catalog No 15-0073
Lot No 21347001-01
Pack Size 100 µg

Type YEATS
Expressed In *E. coli*
Mol. Wgt. 18.9 kDa
Epitope Tag His

Product Description:

YEATS2 YEATS Domain (accession Q9ULM3, amino acids 200-345) expressed in *E. coli* containing a C-terminal 6xHis tag. The YEATS2 protein is part of the ATAC complex, which has histone acetyltransferase activity on histones H3 and H4 [1]. YEATS2 YEATS Domain recognizes and binds H3K27cr [2].

Formulation:

Recombinant 6xHis-tagged protein at 0.33 mg/mL in 50 mM HEPES pH 7.5, 500 mM NaCl, 10% glycerol, 2 mM TCEP.

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:

YEATS2 YEATS Domain is useful for protein binding and screening experiments examining crotonylated protein substrates.

References:

- [1] Wang et al (2008) *J Biol Chem* 283: 33808-15.
- [2] Zhao et al (2016) *Cell Res* 26: 629-632.

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

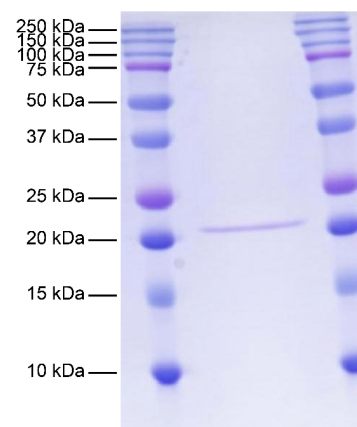


Figure 1: Protein gel data. YEATS2 YEATS Domain (1 µg) was resolved via PAGE gel and stained with Coomassie blue to demonstrate the size and purity of the protein. The migration and molecular weight of the protein standards are indicated.

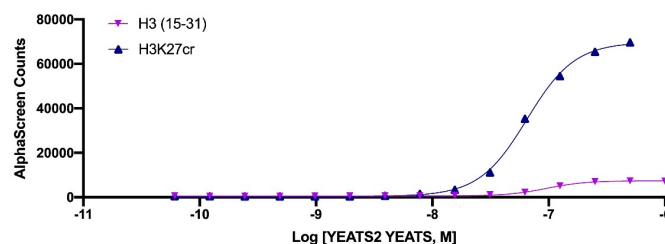


Figure 2: Protein interaction data. YEATS2 YEATS Domain demonstrates binding to H3K27cr peptide and no significant binding to H3(15-31) control peptide when analyzed by AlphaScreen® (PerkinElmer). All peptides assayed at 100 nM.