

GAS41 YEATS Domain, Recombinant Human



EpiCypher®

Catalog No 15-0075
Lot No 22042001-81
Pack Size 100 µg

Type YEATS
Expressed In *E. coli*
Mol. Wgt. 19.5 kDa
Epitope Tag His

Product Description:

GAS41 YEATS Domain (accession O95619, amino acids 15-160) expressed in *E. coli* containing an N-terminal 6xHis tag. The GAS41 protein is part of the NuA4 histone acetyltransferase complex, which is involved with the activation of genes via acetylation of histones H4 and H2A [1]. GAS41 YEATS Domain recognizes and binds acetylated and crotonylated histone H3 [2].

Formulation:

Recombinant 6xHis-tagged protein at 0.90 mg/mL in 20 mM Tris HCl pH 7.5, 500 mM NaCl, 5% glycerol, 1 mM DTT.

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:

GAS41 YEATS Domain is useful for protein binding and screening experiments examining acetylated and crotonylated protein substrates.

References:

- [1] Doyon et al (2004) *Mol Cell Biol* 24: 1884-1896.
[2] Cho et al (2018) *ACS Chem Biol* 13: 2739-2746.

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

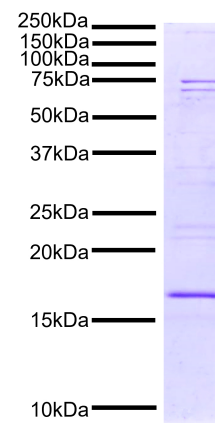


Figure 1: Protein gel data. GAS41 YEATS Domain (1 µg) was resolved via PAGE gel and stained with Coomassie blue to demonstrate the size and purity of the protein. The migration and molecular weight of the protein standards are indicated.

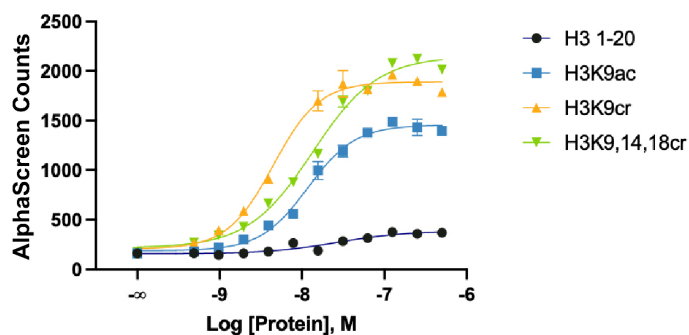


Figure 2: Protein interaction data. GAS41 YEATS Domain demonstrates preferential binding to H3K9cr, followed by H3K9,14,18cr and H3K9ac. There was no significant binding to H3(1-20) control peptide when analyzed by AlphaScreen® (PerkinElmer). All peptides assayed at 100 nM.