

Recombinant YEATS2 (200-340) protein

Catalog No: 81099, 81899

Lot No: 34217001

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 0.9 µg/µl

Source: Human

Buffer Contents: Recombinant YEATS2 (200-340) protein is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, 10% glycerol and 0.5 mM TCEP.

Background: YEATS2 (YEATS Domain Containing 2), is a component of the ATAC complex, a complex with histone acetyltransferase activity on histones H3 and H4. YEATS2 specifically recognizes and binds histone H3 crotonylated at 'Lys-27' (H3K27cr). Crotonylation marks active promoters and enhancers and confers resistance to transcriptional repressors.

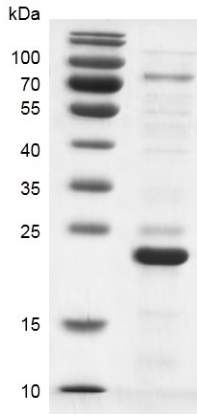
Protein Details: Recombinant human YEATS2 (200-340) that includes amino acids 200-340 (accession number NP_001338299.1) was expressed in *E. coli* cells with an N-terminal 6×His tag and a C-terminal Flag tag. The molecular weight of the protein is 21.5 kDa.

Application Notes: Recombinant YEATS2 (200-340) protein is suitable for use in enzyme kinetics, inhibitor screening, and selectivity profiling.

Binding Assay Conditions: 3 µM H3K27cr (crotonylated Lysine 27) peptide was incubated with different concentrations of YEATS2 (200-340) protein in a 10 µl reaction system containing 50mM HEPES-NaOH pH7.5, 0.1% BSA for 1 hour, then 10 µl FLAG antibody and SA-XL665 mixture (each 1:100 dilution in Binding Buffer) was added to each reaction system and incubated for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

YEATS2 (200-340)



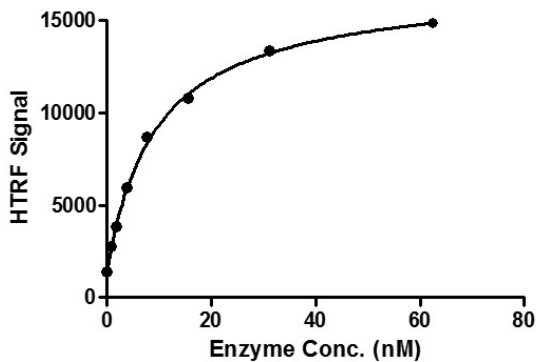
Recombinant YEATS2 (200-340) protein gel

12.5% SDS-PAGE Coomassie staining

MW: 21.5 kDa

Purity: >80%

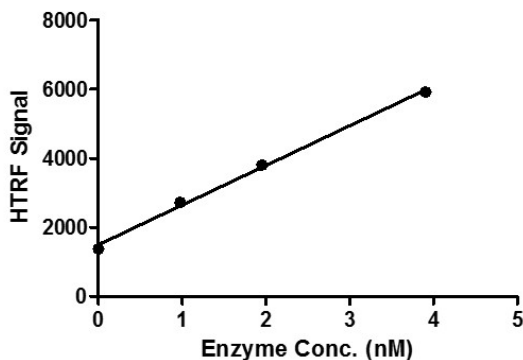
YEATS2 (200-340) Titration



Recombinant YEATS2 (200-340) protein activity assay

H3K27cr (crotonylated Lysine 27) peptide was incubated with different concentrations of YEATS2 (200-340) protein in reaction buffer for 1 hour, then FLAG antibody and SA-XL665 mixture was added to each reaction system for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.

YEATS2 (200-340) Titration



Recombinant YEATS2 (200-340) protein activity assay

H3K27cr (crotonylated Lysine 27) peptide was incubated with different concentrations of YEATS2 (200-340) protein in reaction buffer for 1 hour, then FLAG antibody and SA-XL665 mixture was added to each reaction system for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.