

KAT3B (EP300) Protein

Recombinant human protein expressed in Sf9 cells

Catalog # K312-31BG

Lot # E012-1

Product Description

Recombinant human KAT3B (EP300) (532-1153; contains the catalytic domain) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_001429](#).

Gene Aliases

KAT3B; EP300; p300

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

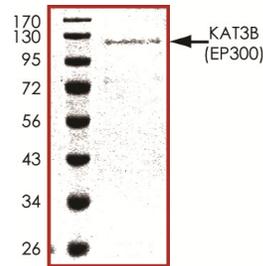
KAT3B (also known as EP300) encodes the adenovirus E1A-associated cellular p300 transcriptional co-activator protein that functions as a histone acetyltransferase. KAT3B regulates transcription via chromatin remodeling and is important in the processes of cell proliferation and differentiation. KAT3B binding is a highly accurate process and can be used for identifying enhancers and their associated activities (1). KAT3B is also useful in the study of the role of tissue-specific enhancers in human biology and diseases on a genome wide scale. Defects in KAT3B gene are a cause of Rubinstein-Taybi syndrome that play a important role in epithelial cancer and act as a classic tumor suppressor gene (2).

References

1. Visel, A. et.al: ChIP-seq accurately predicts tissue-specific activity of enhancers. *Nature* 457: 854-858, 2009.
2. Gayther, S. A. et.al: Mutations truncating the EP300 acetylase in human cancers. *Nature Genet.* 24: 300-303, 2000.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Purity



The purity of KAT3B (EP300) was determined to be **>80%** by densitometry. Approx. MW **120kDa**.

KAT3B (EP300) Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number	K312-31BG
Specific Lot Number	E012-1
Purity	>80%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.