

RBBP4 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # R314-30G

Lot # U1595-3

Product Description

Full-length recombinant human RBBP4 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC075836](#).

Gene Aliases

NURF55, RBAP48

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.

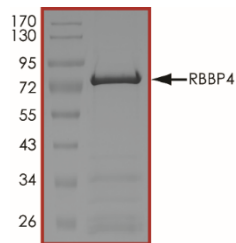
Scientific Background

Histone-binding protein RBBP4 (RBBP4) was initially identified as one of the strongest interactions when a CREB-CREB binding protein (CBP) complex was used as bait to screen a mouse embryo cDNA library in yeast. May target chromatin assembly factors, chromatin remodeling factors and histone deacetylases to their histone substrates in a manner that is regulated by nucleosomal DNA.

References

1. Zhang Q, et al: Histone binding protein RbAp48 interacts with a complex of CREB binding protein and phosphorylated CREB. Mol Cell Biol. 2000; 20(14):4970-8.
2. <http://www.uniprot.org/uniprot/Q09028>

Purity



The purity of RBBP4 was determined to be **>90%** by densitometry. Approx. MW **75 kDa**.

RBBP4 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #	R314-30G
Lot #	U1595-3
Purity	>90%
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.

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

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.