

RBBP5 Protein

Full-length recombinant human protein expressed in E. coli cells

Catalog # R315-30H

Lot # B2093-8

Product Description

Full-length recombinant human RBBP5 was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM_005057](#).

Gene Aliases

RBQ3; SWD1

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 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.

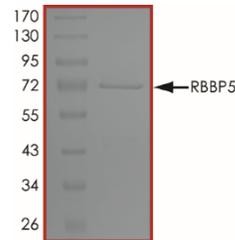
Scientific Background

Retinoblastoma-binding protein 5 (RBBP5) is believed to play a crucial role in the differentiation potential, particularly along the neural lineage in embryonic stem (ES) cells, regulating gene induction and H3 'Lys-4' methylation at key developmental loci, including that mediated by retinoic acid (By similarity). As part of the MLL1/MLL complex, involved in mono-, di- and trimethylation at 'Lys-4' of histone H3. Histone H3 'Lys-4' methylation represents a specific tag for epigenetic transcriptional activation.

References

1. [Thhttp://www.uniprot.org/uniprot/Q15291](http://www.uniprot.org/uniprot/Q15291)
2. Patel A, et al: On the mechanism of multiple lysine methylation by the human mixed lineage leukemia protein-1 (MLL1) core complex. J Biol Chem. 2009; 284(36):24242-56.

Purity



The purity of RBBP5 was determined to be **>95%** by densitometry. Approx. MW **72 kDa**.

RBBP5 Protein

Full-length recombinant human protein expressed in E. coli cells

Catalog #	R315-30H
Lot #	B2093-8
Purity	>95%
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.