

4EBP1 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog # E35-54H

Lot # R145-1

Product Description

Recombinant full-length human 4EBP1 Protein was expressed in Sf9 insect cells using an N-terminal His tag. The gene accession number is [NM_004095](#).

Gene Aliases

BP-1, EIF4EBP1, PHAS-I, MGC4316

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.

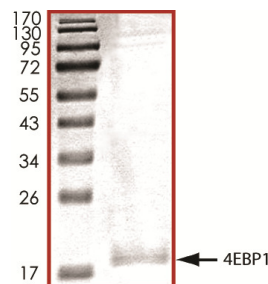
Scientific Background

4EBP1 is a member of a family of translation repressor proteins that directly interact with eukaryotic translation initiation factor 4E (EIF4E). Interaction of 4EBP1 with EIF4E inhibits the multisubunit complex that recruits 40S ribosomal subunits to the 5' end of mRNAs thereby leading to repression of translation. Insulin treatment of adipose cells increases the phosphorylation of 4EBP1 and leads to reduced interaction of 4EBP1 with EIF4E (1). 4EBP1 is expressed in most tissues, with highest levels seen in adipose tissue, pancreas, and skeletal muscle (2).

References

1. Pause, A. et al: Insulin-dependent stimulation of protein synthesis by phosphorylation of a regulator of 5-prime-cap function. *Nature* 371: 762-767, 1994.
2. Tsukiyama-Kohara, K. et al: Tissue distribution, genomic structure, and chromosome mapping of mouse and human eukaryotic initiation factor 4E-binding proteins 1 and 2. *Genomics* 38: 353-363, 1996.

Purity



The purity of 4EBP1 was determined to be **>85%** by densitometry. Approx. MW **18kDa**.

4EBP1 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	E35-54H
Specific Lot Number	R145-1
Purity	>85%
Concentration	0.2 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.