

STUB1 (CHIP) Protein

Full length recombinant protein expressed in E.coli cells

Catalog # C268-30H

Lot # F530-3

Product Description

Recombinant full length human STUB1 (CHIP) was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM_005861](#).

Gene Aliases

CHIP; UBOX1; HSPABP2; NY-CO-7; SDCCAG7

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.

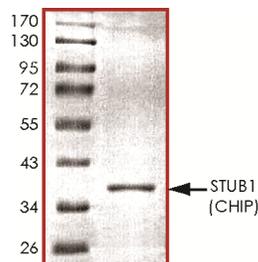
Scientific Background

CHIP(STUB1) or STIP1 homology and U-box containing protein 1 (also known as E3 ubiquitin protein ligase) is a ubiquitin ligase/cochaperone that participates in protein quality control by targeting a broad range of chaperone protein substrates for degradation (1). CHIP acts as a molecular switch for the degradation of misfolded proteins via the proteasome and lysosome pathways (2). CHIP controls the cellular level of BER enzymes and, correspondingly, the efficiency and capacity of BER. CHIP E3 ubiquitin ligases act sequentially in ER membrane and cytosol to monitor the folding status of CFTR and delF508. CHIP possesses intrinsic E3 ubiquitin ligase activity and promotes ubiquitylation.

References

1. Min, J.-N. et.al: CHIP deficiency decreases longevity, with accelerated aging phenotypes accompanied by altered protein quality control. *Molec. Cell. Biol.* 28: 4018-4025, 2008.
2. Shin, Y.et.al: The co-chaperone carboxyl terminus of Hsp70-interacting protein (CHIP) mediates alpha-synuclein degradation decisions between proteasomal and lysosomal pathways. *J. Biol. Chem.* 280: 23727-23734, 2005.

Purity



The purity of STUB1 (CHIP) was determined to be **>95%** by densitometry. Approx. MW **37kDa**.

STUB1 (CHIP) Protein

Full length recombinant protein expressed in E. coli cells

Catalog Number	C268-30H
Specific Lot Number	F530-3
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.