

UBA3 (UBE1C) Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # U203-30G

Lot # F507-1

Product Description

Recombinant full-length human UBA3 (UBE1C) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC022853](#).

Gene Aliases

UBE1C, UBA3, hUBA3

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

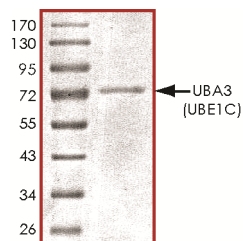
Scientific Background

UBA3 (also known as UBE1C) or ubiquitin-like modifier activating enzyme 3 is a member of the E1 ubiquitin-activating enzyme family. UBA3 associates with AppBp1, an amyloid beta precursor protein binding protein, to form a heterodimer, and then the enzyme complex activates NEDD8, a ubiquitin-like protein, which regulates cell division, signaling and embryogenesis. The mechanism of NEDD8 activation by APPBP1-UBA3 shows a high degree of conservation with ubiquitin activation by UBA1 (1). The NEDD8-activating enzyme, or NAE, composed of NAE1 and UBA3 subunits, is an essential component for the treatment of cancer disease (2).

References

1. Bohnsack, R. N. et.al: Conservation in the mechanism of Nedd8 activation by the human AppBp1-Uba3 heterodimer. J. Biol. Chem. 278: 26823-26830, 2003.
2. Soucy, T. A. et.al: An inhibitor of NEDD8-activating enzyme as a new approach to treat cancer. Nature 458: 732-736, 2009.

Purity



The purity of UBA3 (UBE1C) was determined to be **>95%** by densitometry. Approx. MW **73kDa**.

UBA3 (UBE1C) Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	U203-30G
Specific Lot Number	F507-1
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.