

UBE2U Protein

Full length human recombinant protein expressed in Sf9 cells

Catalog # U237-30G

Lot # O1060-2

Product Description

Recombinant full-length human UBE2U was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_152489](#).

Gene Aliases

RP4-636O23.1

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.

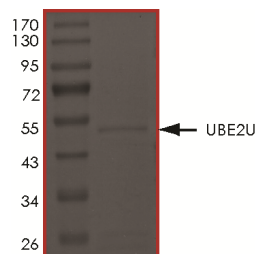
Scientific Background

UBE2U, or Ubiquitin-conjugating enzyme E2 U, belongs to the ubiquitin-conjugating enzyme family. In the global yeast two-hybrid (Y2H) screens to study the specificity of interactions between catalytic domains of the human E2s with RING-type E3s, the analysis showed the UBE2U had a high-quality and physical interactions with a few E3 ligases (1) and (2).

References

1. van Wijk SJ, et.al: A comprehensive framework of E2-RING E3 interactions of the human ubiquitin-proteasome system. Mol Syst Biol. 5:295. doi: 10.1038, 2009.
2. Markson G, et.al: Analysis of the human E2 ubiquitin conjugating enzyme protein interaction network. Genome Res. 19 (10):1905-11. doi: 10.1101, 2009.

Purity



The purity of UBE2U was determined to be **>80%** by densitometry. Approx. MW **52kDa**.

UBE2U Protein

Full length human recombinant protein expressed in Sf9 cells

Catalog Number	U237-30G
Specific Lot Number	O1060-2
Purity	>80%
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.