

UBE2Z (USE1) Protein

Recombinant protein expressed in E.coli cells

Catalog # U241-31H

Lot # J689-4

Product Description

Recombinant human UBE2Z (USE1) (109-end) was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM_023079](#).

Gene Aliases

HOYS7; USE1

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.

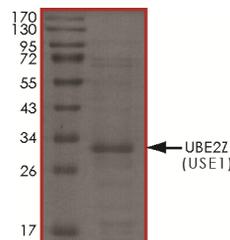
Scientific Background

UBE2Z(USE1) encodes the ubiquitin conjugating enzyme (E2) which is involved in Ubiquitination, the covalent attachment of the protein ubiquitin (Ub) to other cellular proteins (1). UBE2Z plays an important role in the ubiquitin process and has been implicated in a number of important physiological processes including apoptosis. UBE2Z is specifically charged with ubiquitin by UBA6 (2). UBE2Z is widely expressed in human tissues and expression is especially high in placenta, pancreas, spleen and testis.

References

1. Gu, X. et al: Cloning and characterization of a gene encoding the human putative ubiquitin conjugating enzyme E2Z (UBE2Z). Mol Biol Rep. 2007 Sep;34(3):183-8.
2. Jin, J. et.al: Dual E1 activation systems for ubiquitin differentially regulate E2 enzyme charging. Nature 447: 1135-1138, 2007.

Purity



The purity of UBE2Z (USE1) was determined to be **>80%** by densitometry. Approx. MW **30 kDa**.

UBE2Z (USE1) Protein

Recombinant protein expressed in E. coli cells

Catalog Number	U241-31H
Specific Lot Number	J689-4
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.