

UBE2D2 (UBC4) Protein

Full length recombinant protein expressed in E.coli cells

Catalog # U214-30H

Lot # F698-2

Product Description

Recombinant full length *Xenopus* UBE2D2 (UBC4) was expressed in E. coli cells using a N-terminal His tag. The gene accession number is [NM_001093036](#).

Gene Aliases

UBC4; PUBC1; UBC4/5; UBCH5B; E2(17)KB2

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.

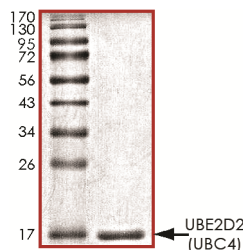
Scientific Background

UBE2D2(UBC4) Protein or ubiquitin-conjugating enzyme E2D 2 is a member of the E2 ubiquitin-conjugating enzyme family which functions in the ubiquitination of the tumor-suppressor protein p53, which is induced by an E3 ubiquitin-protein ligase. UBE2D2 specifically ubiquitinates E6AP and in vivo inhibition of UBE2D2 can lead to inhibition of E6-stimulated p53 degradation (1). UBE2D2 could conjugate ubiquitin to target proteins in an E6AP-dependent manner (2).

References

1. Rolfe, M.et.al: Reconstitution of p53-ubiquitylation reactions from purified components: the role of human ubiquitin-conjugating enzyme UBC4 and E6-associated protein (E6AP). Proc. Nat. Acad. Sci. 92: 3264-3268, 1995.
2. Jensen, J. P.et.al: Identification of a family of closely related human ubiquitin conjugating enzymes. J. Biol. Chem. 270: 30408-30414, 1995.

Purity



The purity of UBE2D2 (UBC4) was determined to be **>95%** by densitometry. Approx. MW **17 kDa**.

UBE2D2 (UBC4) Protein

Full length recombinant protein expressed in E. coli cells

Catalog Number	U214-30H
Specific Lot Number	F698-2
Purity	>95%
Concentration	0.1µg/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.