

Recombinant UBE2D3 protein

Catalog No: 31230

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 0.5 µg/µl

Source: Human

Buffer Contents: 10 µg of recombinant protein supplied at a concentration of 0.5 µg/µl in 20 mM Tris-HCl, pH 7.3, 100 mM KCl, 0.2 mM EDTA and 20% glycerol.

Background: UBE2D3, ubiquitin-conjugating enzyme E2D 3, (also known as UBCH5C, UBC4/5, MGC5416, MGC43926, or E2(17)KB3) belongs to the E2 ubiquitin-conjugating enzyme family that covalently attaches ubiquitin to other proteins and marks them for degradation. Ubiquitination involves three classes of enzymes: E1s, ubiquitin-activating enzymes, E2s, ubiquitin-conjugating enzymes and E3s, ubiquitin-protein ligases. UBE2D2 functions in the E6/E6-AP-induced ubiquitination of the tumor suppressor protein p53.

Protein Details: Recombinant UBE2D3 protein was expressed in *E. coli* as the full-length human protein (accession number AAA91461) with an N-terminal poly-histidine tag. The molecular weight of the protein is ~18.7 kDa.

Application Notes: Recombinant UBE2D3 protein is suitable for use in ubiquitination assays. A recommended starting point for the ubiquitination assay is 200-500 ng recombinant protein per reaction.

Storage and Guarantee: This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.