

## Recombinant UBE2M protein

---

**Catalog No:** 31235

**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:** UBE2M, ubiquitin-conjugating enzyme E2M (also known as UBC12, ubiquitin-conjugating enzyme E2M (UBC12 homolog, yeast) or Nedd8-conjugating enzyme Ubc12) is an E2 ubiquitin-conjugating enzyme that is linked with NEDD8 and can be conjugated to cellular proteins. It accepts the ubiquitin-like protein NEDD8 from the UBA3-NAE1 E1 complex and catalyzes covalent attachment to other proteins.

**Protein Details:** Recombinant UBE2M protein was expressed in *E. coli* as the full-length human protein (accession number AAC26141) with an N-terminal poly-histidine tag. The molecular weight of the protein is ~23 kDa.

**Application Notes:** Recombinant UBE2M protein is suitable for use in neddylation assays. A recommended starting point for the neddylation 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.