

## UBE2D3 recombinant protein

**Other names:** UBCH5C, UBC5C, PR02116, UBC4/5

**Cat. No.** C23010112

**Source:** E. coli

**Lot #:** 001

**Size:** 25 µg/ 25 µl

**Concentration:** 1 µg/µl

**Specificity:** Human

**Purity:** Purified using FPLC, >98% purity as determined by SDS-PAGE.

**Storage buffer:** 20 mM Tris-Cl pH 7.6, 0.15 M NaCl, 10% glycerol, and 1 mM DTT.

**Storage:** Store at -80°C; guaranteed stable for 2 years from date of receipt when stored properly.

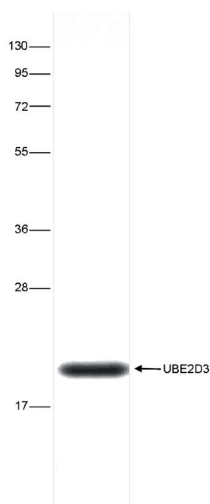
**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**Description:** Full length recombinant human N-terminal His tagged UBE2D3 (Ubiquitin-Conjugating Enzyme E2 3) protein, produced in E. coli.

### Protein description

UBE2D3 (UniProt/Swiss-Prot entry P61077) is a member of the E2 ubiquitin-conjugating enzyme family. It accepts ubiquitin from the E1 complex and catalyzes its covalent attachment to target proteins.

### Quality control



**Figure 1.**

SDS page of the UBE2D3 recombinant protein. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.

#### Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing (Ougrée) - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

#### Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

**Last update:** August 28, 2015