

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 E2D 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

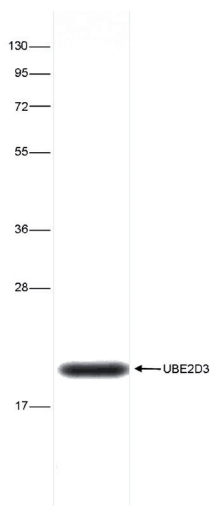


Figure 1.

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