

Ubiquitin recombinant protein

Cat. No. C23010113

Source: E. coli

Lot #: 001

Size: 25 µg/ 50 µl

Concentration: 0.5 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 Ubiquitin, produced in E. coli and biotinylated in vitro.

Protein description

Covalent binding of ubiquitin to other proteins is one of the most frequent posttranslational modifications. This covalent attachment can be either as a monomer (monoubiquitin), a polymer linked via different Lys residues of the ubiquitin (polyubiquitin chains) or a linear polymer linked via the initiator Met of the ubiquitin (linear polyubiquitin chains). Conjugation of ubiquitin to its target protein can lead to various effects within a cell, including protein degradation, DNA repair, cell cycle regulation, kinase modification, endocytosis, and regulation of other cell signaling pathways.

Quality control

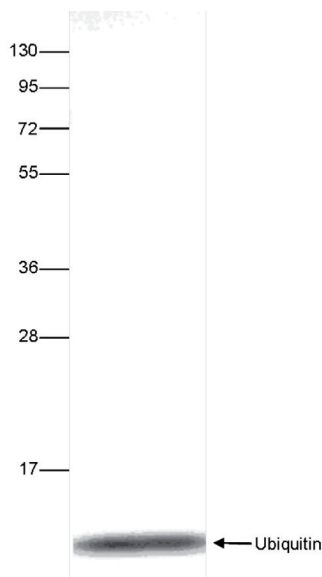


Figure 1.

SDS page of the biotinylated Ubiquitin 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 31, 2015