

Recombinant Ubiquitin (1-76) protein, His-Tag

Catalog No: 82018, 82618

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Source: Human

Buffer Contents: Recombinant Ubiquitin (1-76) protein is supplied in 25 mM Tris 8.0, 300mM NaCl, 20% glycerol, 0.5 mM TCEP.

Background: Ubiquitin exists either covalently attached to another protein, or free (unanchored). When covalently bound, it is conjugated to target proteins via an isopeptide bond 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). Polyubiquitin chains, when attached to a target protein, have different functions depending on the Lys residue of the ubiquitin that is linked. Linear polymer chains formed via attachment by the initiator Met lead to cell signaling. **Ubiquitin** is usually conjugated to Lys residues of target proteins, however, in rare cases, conjugation to Cys or Ser residues has been observed. When polyubiquitin is free (unanchored-polyubiquitin), it also has distinct roles, such as in activation of protein kinases, and in signaling.

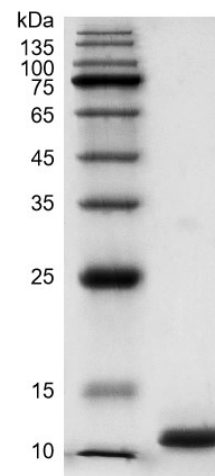
Protein Details: Recombinant Ubiquitin (1-76) protein that includes amino acids 1-76 of human UBC protein (accession number NP_066289.3) was expressed in *E. coli* and contains an N-terminal His tag with a molecular weight of 10.73 kDa. The purity of the protein is ≥ 95% by SDS-PAGE

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for this product is shown.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. 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.

Ubiquitin(1-76)



Recombinant Ubiquitin (1-76) protein
12.5% SDS-PAGE Coomassie staining
MW: 10.73 kDa
Purity: >95%