

## Recombinant Ubiquitin protein

---

**Catalog No:** 31231

**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:** Ubiquitin is a small, highly conserved regulatory protein that exists in all eukaryotic cells. Ubiquitin is covalently attached to a range of target proteins in order to mark these proteins for proteasomal degradation. Ubiquitination involves three classes of enzymes: E1s, ubiquitin-activating enzymes, E2s, ubiquitin-conjugating enzymes and E3s, ubiquitin-protein ligases.

**Protein Details:** Recombinant Ubiquitin protein was expressed in *E. coli* as amino acids 1-76 of human protein (accession number AAA72608) with an N-terminal poly-histidine tag. The molecular weight of the protein is ~10.6 kDa.

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