

## Recombinant Ubiquitin G76A protein

---

**Catalog No:** 31232

**Source:** Human

**Quantity:** 10 µg

**Expressed In:** *E. coli*

**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 G76A protein was expressed in *E. coli* as amino acids 1-76 of human protein (accession number AAA72608) with an N-terminal poly-histidine tag. Recombinant ubiquitin G76A protein contains a substitution of a glycine residue to alanine at amino acid 76. The molecular weight of the protein is ~10.6 kDa.

### Protein Applications:

Recombinant Ubiquitin G76A protein is suitable for use in ubiquitination assays as a negative control. A recommended starting point for the ubiquitination assay is 200-500 ng recombinant protein per reaction.

### Storage:

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

### Guarantee:

This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.