

Recombinant NEDD8 G76A protein

Catalog No: 31237

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:

NEDD8, neural precursor cell expressed, developmentally down-regulated 8 (also known as **NEDD-8**, **FLJ43224**, **MGC104393**, **MGC125896**, or **MGC125897**) is a ubiquitin-like protein (Ubl) that regulates cell division, signaling and embryogenesis. NEDD8 is activated by the heterodimeric ubiquitin activating enzyme E1 complex, UBA3-NAE1 E1, and the E2 enzyme UBE2M catalyzes the covalent attachment to target proteins in the presence of the appropriate E3 ligase. There are several proteases capable of removing NEDD8 from its protein conjugates.

Protein Details:

Recombinant NEDD8 G76A protein was expressed in *E. coli* as amino acids 1-76 of human protein (accession number NP_006147) with an N-terminal poly-histidine tag. Recombinant NEDD8 G76A protein contains a substitution of a glycine residue to alanine at amino acid 76. The molecular weight of the protein is ~10.5 kDa.

Protein Applications:

Recombinant NEDD8 G76A protein is suitable for use in neddylation assays as a negative control. A recommended starting point for the neddylation 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.