

Recombinant NEDD8 protein

Catalog No: 31236

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: 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 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. The molecular weight of the protein is ~10.5 kDa.

Application Notes: Recombinant NEDD8 protein is suitable for use in neddylation assays. A recommended starting point for the neddylation 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.