

Recombinant NAE1 protein

Catalog No: 31233

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: NAE1, NEDD8 activating enzyme E1 subunit 1 (also known as APP-BP1, APPBP1, or amyloid beta precursor protein-binding protein 1) is a subunit of the heterodimeric ubiquitin activating enzyme E1, which is specific for the Rub family of ubiquitin-like proteins (Ubls). NAE1 associates with UBA3 to form the UBA3-NAE1 E1 enzyme complex that activates the ubiquitin-like protein NEDD8. NEDD8 regulates cell division, signaling and embryogenesis.

Protein Details: Recombinant NAE1 protein was expressed in *E. coli* as the full-length human protein (accession number AAC50477) with an N-terminal poly-histidine tag. The molecular weight of the protein is ~62.3 kDa.

Application Notes: Recombinant NAE1 protein is suitable for use in neddylation assays. A recommended starting point for the neddylation assay is 100 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.