

Recombinant CDKN1B protein

Catalog No: 81390, 81690

Expressed In: *E. coli*

Quantity: 50, 1000 µg

Concentration: 0.5 µg/µl

Source: Human

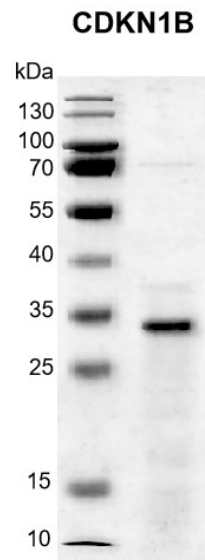
Buffer Contents: Recombinant CDKN1B protein is supplied in 25 mM Tris pH 8.0, 300 mM NaCl, 10% glycerol, 0.5 mM TCEP.

Background: CDKN1B is a cyclin-dependent kinase inhibitor, also called cyclin-dependent kinase inhibitor p27 or p27Kip1, important regulator of cell cycle progression. Inhibits the kinase activity of CDK2 bound to cyclin A, but has little inhibitory activity on CDK2 bound to SPDYA. Involved in G1 arrest. Potent inhibitor of cyclin E- and cyclin A-CDK2 complexes. Forms a complex with cyclin type D-CDK4 complexes and is involved in the assembly, stability, and modulation of CCND1-CDK4 complex activation. Acts either as an inhibitor or an activator of cyclin type D-CDK4 complexes depending on its phosphorylation state and/or stoichiometry.

Protein Details: Recombinant CDKN1B protein was expressed in *E. coli* as the full length protein (accession number NP_004055.1) with an N-terminal 6×His tag. The molecular weight of the protein is 24.01 kDa

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of arrival.



Recombinant CDKN1B protein
12.5% SDS-PAGE Coomassie staining

M.W.: 24.01 kDa

Purity: ≥ 80%