

Recombinant NAP1L1 protein

Catalog No: 31508, 31912

Expressed In: *E. coli*

Quantity: 50, 1000 µg

Concentration: 5 µg/µl

Source: Human

Buffer Contents: Full length recombinant NAP1L1 protein expressed in *E. coli* cells and supplied at a concentration of 5 µg/µl in 20 mM Tris pH 8.0, 100 mM KCl, 20% Glycerol, 0.2 mM EDTA.

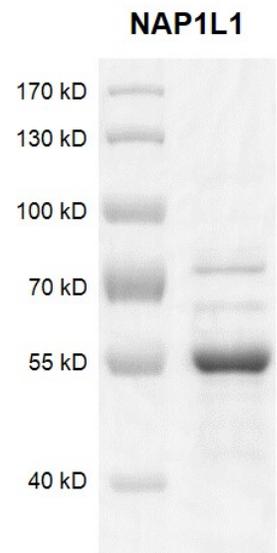
Background: **NAP1L1** (Nucleosome assembly protein 1-like 1; NAP1-Related Protein, NRP) is a member of the NAP family of histone chaperones. Histone chaperones are involved in chromatin assembly following DNA replication and may play a role in modulating chromatin formation and contribute to the regulation of cell proliferation.

Protein Details: Recombinant NAP1L1 protein (accession number NP_004528.1) was expressed in *E. coli* cells and contains an N-terminal His tag with a molecular weight of 47.5 kDa. The recombinant protein is >80% pure by SDS-PAGE.

Application Notes: Recombinant NAP1L1 is suitable for in vitro chromatin assembly and protein-protein interaction.

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 receipt.

This product is for research use only and is not for use in diagnostic procedures.



Recombinant NAP1L1 protein gel. NAP1L1 protein was run on an 8% SDS-PAGE gel and stained with Coomassie blue.