

Recombinant CENPB protein

Catalog No: 81201, 81901

Lot No: 24718001

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 0.3 µg/µl

Source: Human

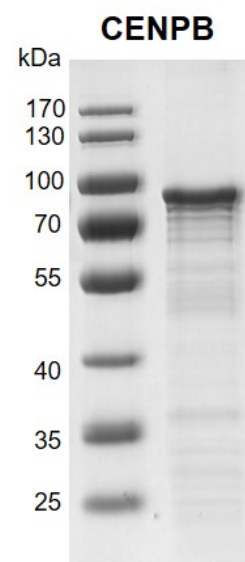
Buffer Contents: Recombinant CENPB protein is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, 10% glycerol, and 0.5 mM TCEP.

Background: CENPB (Centromere Protein B), also called as CENP-B, is a highly conserved protein that facilitates centromere formation. It is a DNA-binding protein that is derived from transposases of the pogo DNA transposon family. It contains a helix-loop-helix DNA binding motif at the N-terminus, and a dimerization domain at the C-terminus. CENPB can interact with centromeric heterochromatin in chromosomes and binds to a specific 17 bp subset of alphoid satellite DNA, called the CENP-B box. It may organize arrays of centromere satellite DNA into a higher-order structure which then directs centromere formation and kinetochore assembly in mammalian chromosomes. CENPB is proposed to play an important role in the assembly of specific centromere structures in interphase nuclei and on mitotic chromosomes. It is also considered a major centromere autoantigen recognized by sera from patients with anti-centromere antibodies.

Protein Details: Recombinant human CENPB protein was expressed in *E. coli* cells as the full length protein (accession number NP_001801) with an N-terminal 6xHis -Tag. The molecular weight of the protein is 68.7 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 this product is shown.

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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



Recombinant CENPB protein gel

10% SDS-PAGE Coomassie staining
MW: 68.7 kDa
Purity: >85%