

Recombinant RANBP2 (2553-2838) protein

Catalog No: 82101, 82701

Quantity: 100, 1000 µg

Expressed In: *E. coli*

Source: Human

Buffer Contents: Recombinant RANBP2 (2553-2838) protein is supplied in 25 mM Tris pH 8.0, 300 mM NaCl, 20% glycerol, 0.5 mM TCEP.

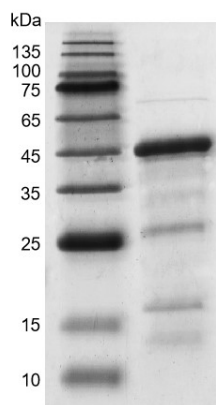
Background: **RANBP2 (E3 SUMO-protein ligase RanBP2)** is a E3 SUMO-protein ligase which facilitates SUMO1 and SUMO2 conjugation by UBE2I. Involved in transport factor (Ran-GTP, karyopherin)-mediated protein import via the F-G repeat-containing domain which acts as a docking site for substrates. Could also have isomerase or chaperone activity and may bind RNA or DNA. Component of the nuclear export pathway. Specific docking site for the nuclear export factor exportin-1.

Protein Details: Recombinant RANBP2 (2553-2838) protein that includes amino acids 2553-2838 of human RANBP2 protein (accession number NP_006258.3) was expressed in *E. coli* and contains an N-terminal His tag with a molecular weight of 34.05 kDa. The purity of the protein is ≥ 90% by SDS-PAGE.

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

RANBP2 (2553-2838)



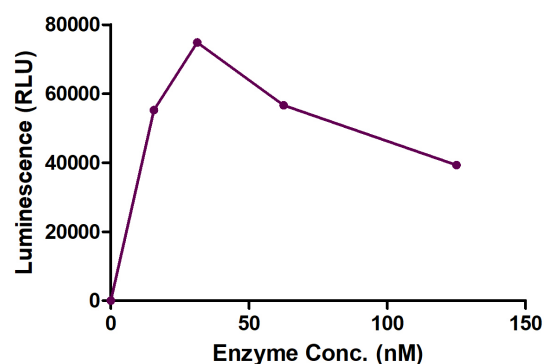
Recombinant RANBP2 (2553-2838) protein gel

12.5% SDS -PAGE Coomassie staining

MW: 34.05 kDa

Purity: $\geq 80\%$

RANBP2 (2553-2838) Titration



ELISA for RANBP2 (2553-2838) activity

ELISA plate was coated with 50 ng p53 protein in each well. Then 100 nM SAE1/SAE2 Complex, 200 nM UBE2I, 5 μ M SUMO2 -biotin, 100 μ M ATP and different concentrations of RANBP2 protein in 50 μ l reaction system containing 40 mM Tris-HCl pH 7.4, 20 mM MgCl₂, 1 mM DTT, 0.1 mg/ml BSA were added in each well and incubated at 37°C for 1 hour. Strep-HRP was used to detect reaction products.