

Recombinant UBE2L6 protein

Catalog No: 81400, 81500

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Source: Human

Buffer Contents: Recombinant UBE2L6 protein is supplied in 25 mM Tris 8.0, 300mM NaCl, 20% glycerol, and 0.5 mM TCEP.

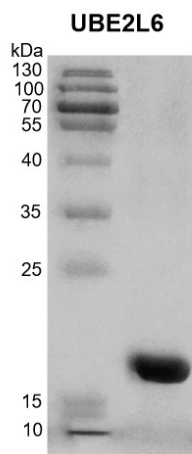
Background: UBE2L6 (Ubiquitin/ISG15-conjugating enzyme E2 L6) also known as RIG-B and UBCH8. UBE2L6 can catalyze the covalent attachment of ubiquitin or ISG15 to other proteins. Functions in the E6/E6-AP-induced ubiquitination of p53/TP53. Promotes ubiquitination and subsequent proteasomal degradation of FLT3.

Protein Details: Recombinant UBE2L6 protein that includes full length of human UBE2L6 protein (accession number NP_004214.1) was expressed in *E. coli* and contains an N-terminal His tag with a molecular weight of 19.93 kDa. The purity of the protein is ≥ 95% 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 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 guaranteed for 6 months from date of receipt.

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

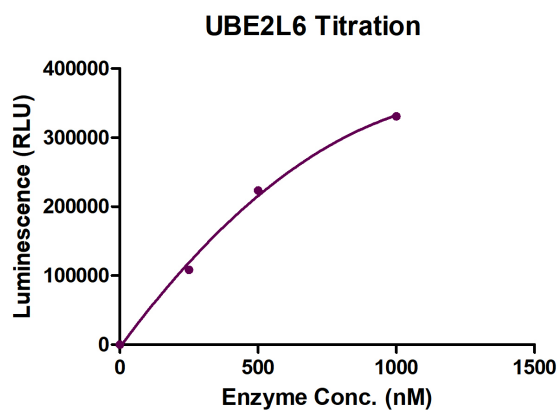


Recombinant UBE2L6 protein gel.

Recombinant UBE2L6 was run on a 12.5% SDS-PAGE gel and stained with Coomassie Blue.

MW: 19.93 kDa

Purity: >95%



AMP-Glo assay for UBE2L6 activity

7.9 μ M ubiquitin, 63 nM UBA1 and 25 μ M ATP were incubated with different concentrations of UBE2L6 in 10 μ l reaction system containing 40 mM Tris-HCl pH 7.4, 20 mM MgCl₂, 0.5 mM DTT, 0.1 mg/ml BSA at 37°C for 1 hour. 10 μ l of AMP-Glo Reagent I was added to the reaction and incubated for 1 hour at room temperature. Then 20 μ l of AMP-Glo Detection Solution was added and luminescence were read after another 30 min incubation.