

Recombinant UBE2L3 protein

Catalog No: 81499, 81599

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Source: Human

Buffer Contents: Recombinant UBE2L3 protein is supplied in 25 mM Tris 8.0, 300mM NaCl, 20% glycerol, 0.5 mM TCEP.

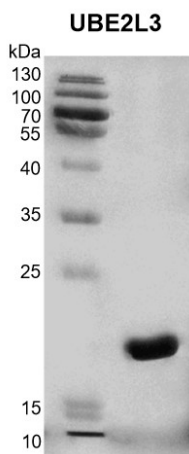
Background: UBE2L3 (Ubiquitin-conjugating enzyme E2 L3) also known as E2-F1, L-UBC, UBCH7 and UbcM4. **UBE2L3** that specifically acts with HECT-type and RBR family E3 ubiquitin-protein ligases. It has activity with the RBR family E3 enzymes, such as PRKN and ARIH1, that function like RING-HECT hybrids. Accepts ubiquitin from the E1 complex and catalyzes its covalent attachment to other proteins. In vitro catalyzes 'Lys-11'-linked polyubiquitination. Involved in the selective degradation of short-lived and abnormal proteins. Down-regulated during the S-phase it is involved in progression through the cell cycle. Regulates nuclear hormone receptors transcriptional activity. May play a role in myelopoiesis.

Protein Details: Recombinant UBE2L3 protein that includes full length of human UBE2L3 protein (accession number NP_003338.1) was expressed in *E. coli* and contains an N-terminal His tag with a molecular weight of 20.02 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 for 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.

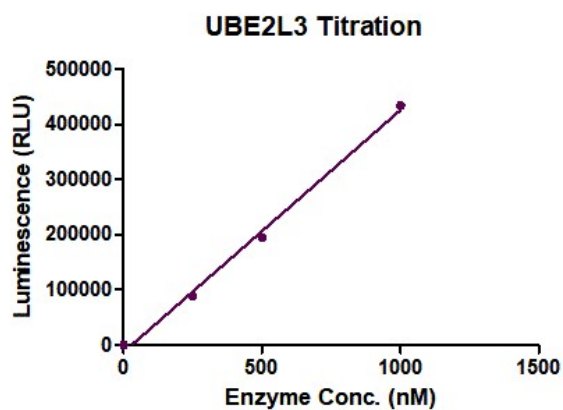


Recombinant UBEL3 protein gel.

12.5% SDS-PAGE Coomassie staining

MW: 20.02 kDa

Purity: $\geq 95\%$



AMP-Glo assay for UBE2L3 activity

7.9 μM ubiquitin, 63 nM UBA1 and 25 μM ATP were incubated with different concentrations of UBE2L3 in 10 μl reaction system containing 40 mM Tris-HCl pH 7.4, 20 mM MgCl_2 , 0.5 mM DTT, 0.1 mg/ml BSA at 37°C for 1 hour. 10 μl of AMP-Glo Reagent I was added to the reaction and incubated for 1 hour at room temperature. Then 20 μl of AMP-Glo Detection Solution was added and luminescence was read after another 30 min incubation.