

Recombinant UBE2D4 protein

Catalog No: 82004, 82604

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Source: Human

Buffer Contents: Recombinant UBE2D4 protein is supplied in 25 mM Tris 7.4, 300mM NaCl, 20% glycerol, 0.5 mM TCEP.

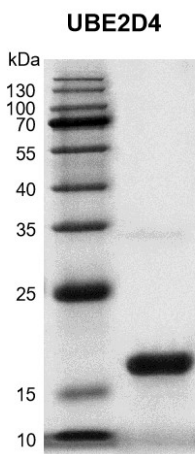
Background: UBE2D4 (Ubiquitin-conjugating enzyme E2 D4) also known as HBUCE1. UBE2D4 is a member of the E2 ubiquitin-conjugating enzyme family which is required for post-replicative DNA damage repair and plays an important role in various cellular processes. UBE2D4 can accept ubiquitin from the E1 complex and catalyzes its covalent attachment to other proteins. In vitro able to promote polyubiquitination using all 7 ubiquitin Lys residues but may prefer 'Lys-11' and 'Lys-48'-linked polyubiquitination.

Protein Details: Recombinant UBE2D4 protein that includes full length of human UBE2D4 protein (accession number NP_057067.1) was expressed in *E. coli* and contains an N-terminal His tag with a molecular weight of 18.81 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.

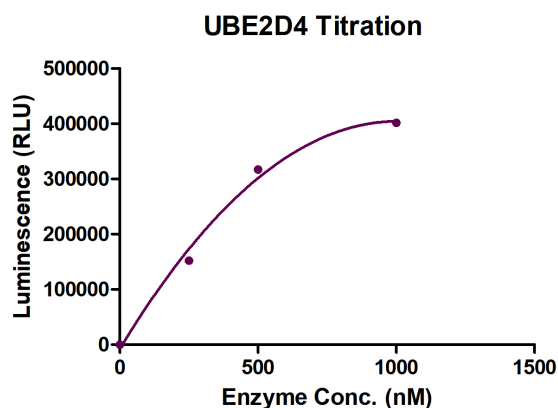


Recombinant UBE2D4 protein gel.

UBE2D4 protein was run on a 12.5% SDS-PAGE gel and stained with Coomassie Blue.

MW: 18.81 kDa

Purity: >95%



AMP-Glo assay for UBE2D4 activity

7.9 μ M ubiquitin, 63 nM UBA1 and 25 μ M ATP were incubated with different concentrations of UBE2D4 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.