

## Recombinant UBE2C protein

**Catalog No:** 82010, 82610

**Expressed In:** *E. coli*

**Quantity:** 100, 1000 µg

**Source:** Human

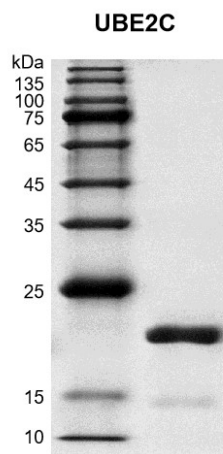
**Buffer Contents:** Recombinant UBE2C protein is supplied in 25 mM Tris 8.0, 300mM NaCl, 20% glycerol, 0.5 mM TCEP.

**Background:** UBE2C (Ubiquitin-conjugating enzyme E2 C) also known as dJ447F3.2 and UBCH10. UBE2C 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. UBE2C can accept ubiquitin from the E1 complex and catalyzes its covalent attachment to other proteins. In vitro catalyzes 'Lys-11'- and 'Lys-48'-linked polyubiquitination. Acts as an essential factor of the anaphase promoting complex/cyclosome (APC/C), a cell cycle-regulated ubiquitin ligase that controls progression through mitosis. Acts by initiating 'Lys-11'-linked polyubiquitin chains on APC/C substrates, leading to the degradation of APC/C substrates by the proteasome and promoting mitotic exit.

**Protein Details:** Recombinant UBE2C protein that includes full length of human UBE2C protein (accession number NP\_008950.1) was expressed in *E. coli* and contains an N-terminal His tag with a molecular weight of 21.82 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 this product is shown.

**Storage and Guarantee:**

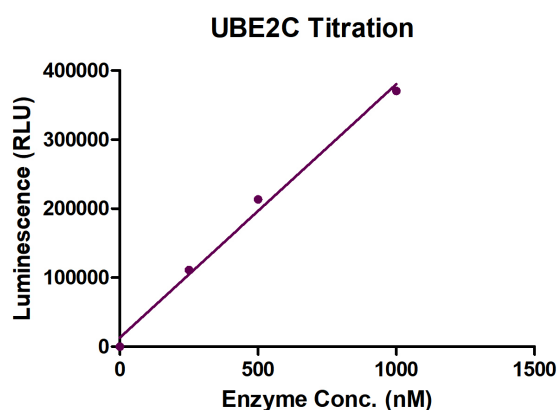


#### Recombinant UBE2C protein gel.

UBE2C protein was run on a 12.5% SDS-PAGE gel and stained with Coomassie Blue.

MW: 21.82 kDa

Purity: >90%



#### AMP-Glo assay for UBE2C activity

7.9  $\mu$ M ubiquitin, 63 nM UBA1 and 25  $\mu$ M ATP were incubated with different concentrations of UBE2C in 10  $\mu$ l reaction system containing 40 mM Tris-HCl pH 7.4, 20 mM MgCl<sub>2</sub>, 0.5 mM DTT, 0.1 mg/ml BSA at 37°C for 1 hour. 10  $\mu$ l of AMP-Glo Reagent I was added to the reaction and incubated for 1 hour at room temperature. Then 20  $\mu$ l of AMP-Glo Detection Solution was added and luminescence was read after another 30 min incubation.