

UBE2C, Active

Recombinant full-length human proteins expressed in *E. coli* cells

Catalog # U212-380H

Lot # V2408-8

Product Description

Recombinant full-length human UBE2C was expressed in *E. coli* cells using an N-terminal His tag. The UBE2C gene accession number is [NM_007019](#).

Gene Aliases

dJ447F3.2; UBCH10

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

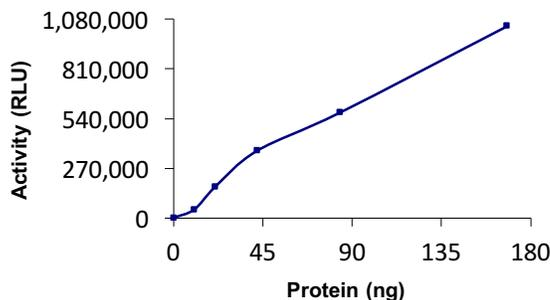
Scientific Background

UBE2C or ubiquitin-conjugating enzyme E2C is a member of the E2 ubiquitin-conjugating enzyme family which is required for the destruction of mitotic cyclins and for cell cycle progression. UBE2C acted as a dominant-negative inhibitor which blocks the destruction of mitotic cyclins A and B and the onset of anaphase in frog embryos and mammalian cells. High UBE2C mRNA expression is associated with poor disease-free survival and overall survival. High tumor grade, as well as high Ki67 protein expression, was more frequent in the high-expression group of UBE2C.

References

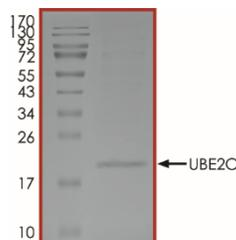
1. Mondal, S. et al: A bioluminescent assay for monitoring conjugation of ubiquitin and ubiquitin-like proteins. *Anal. Biochem.* 510: 41-51, 2016
2. Rape, M. et al: Autonomous regulation of the anaphase-promoting complex couples mitosis to S-phase entry. *Nature* 432: 588-595, 2004
3. Townsley, F. M. et al: Dominant-negative cyclin-selective ubiquitin carrier protein E2-C/Ubch10 blocks cells in metaphase. *Proc. Nat. Acad. Sci.* 94: 2362-2367, 1997

Specific Activity



The specific activity of UBE2C was determined to be **16 nmol/min/mg** as per activity assay protocol.

Purity



The purity of UBE2C was determined to be **>95%** by densitometry, approx. MW **21 kDa**.

UBE2C, Active

Recombinant full-length human protein expressed in *E. coli* cells

Catalog #	U212-380H
Specific Activity	16 nmol/min/mg
Lot #	V2408-8
Purity	>95%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.

Activity Assay Protocol

Reaction Components

Active Ubiquitinating Enzymes

Active UBE2C (Catalog #: U212-380H), UBA1 (Catalog #: U201-380G) and BIRC3 (Catalog #: B280-380G) diluted with Ubiquitination Buffer and assayed as outlined in sample activity plot. (Note: these are suggested working dilutions and it is recommended that the researcher perform a serial dilution of Active UBE2C for optimal results).

Ubiquitination Buffer

Buffer components: 40mM Tris (pH7.5), 20mM MgCl₂, 0.1mg/ml BSA. Add 0.5mM DTT prior to use.

AMP-Glo™ Assay (Promega, Catalog #: V5011)

AMP, 10 mM
Ultra Pure ATP, 10mM
AMP-Glo™ Reagent I
AMP-Glo™ Reagent II
Kinase-Glo™ One Solution

Substrate (Catalog #: U06-54N)

Wild-type ubiquitin protein diluted with Ubiquitination Buffer to a working stock of 170ng/μl (20μM).

Assay Protocol

The UBE2C assay is performed using the AMP-Glo™ Assay kit (Promega), by detecting the amount of the universal AMP generated. Ubiquitin conjugation is proportional to the generated AMP, and the presence of all components of the Ub conjugation machinery (Ub, E1, E2, and E3) is required for maximal activity of the system.

- Step 1.** Thaw the active UBE2C, UBA1, BIRC3 and ubiquitin on ice, and all AMP-Glo™ components except AMP-Glo™ Reagent II at room temperature. Keep AMP-Glo™ Reagent II on ice.
- Step 2.** Prepare the following working solutions with Ubiquitination Buffer:
 - o 2X Reaction Cocktail: 170ng/μl ubiquitin + 15ng/μl UBA1 + 40ng/μl BIRC3 + 50μM ATP
 - o 2X final concentration of Active UBE2C
- Step 3.** In a half-area white 96-well plate, add the following components to bring the initial reaction volume to 10 μl:
 - Component 1.** 5 μl of 2X Reaction Cocktail
 - Component 2.** 5 μl of 2X Active UBE2C

Note: A blank control can be set up as outlined above by replacing the enzyme working solution with an equal volume of Ubiquitination Buffer.
- Step 4.** Briefly centrifuge the plate to ensure reagents are fully mixed and at the bottom of the wells. Seal the plate with a plate seal and incubate at 37°C for 60 minutes
- Step 5.** Equilibrate plate to room temperature. Add 10 μl of AMP-Glo™ Reagent I to all wells, mix by shaking for 1-2 minutes. Incubate the plate at room temperature for 60 minutes.
- Step 6.** Prepare AMP Detection Solution by adding AMP-Glo™ Reagent II to Kinase-Glo™ One Solution at a 1:100 volume ratio. Add 20 μl of the Detection Solution to all wells. Mix for 1-2 minutes and incubate at room temperature for 30 minutes
- Step 7.** Read the plate using the KinaseGlo Luminescence Protocol on a GloMax plate reader (Promega; Cat# E7031)
- Step 8.** Using the AMP standard curve, determine the concentration of AMP produced (μM) and calculate the enzyme specific activity as outlined below. For a detailed protocol of how to determine AMP amount from RLUs, see AMP-Glo™ Assay protocol at Promega's website: www.promega.com/protocols

Enzyme Specific Activity (SA) (nmol/min/mg)

$$= \frac{[AMP](\mu M) \times \text{Reaction Volume}(\mu l)}{\text{Reaction Time (min)} \times \text{Enzyme Amount (mg)}} \times 10^{-3}$$

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.