

Data Sheet

UBE2C, His-Tag

Human, Recombinant, with N-terminal His-Tag
Catalog #: 79369
Lot #: 171110 **Conc.:** 0.1 mg/ml

Formulated in: 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, and 25% glycerol.

Stability: At least 6 months at -70°C.
 Avoid freeze/thaw cycles.

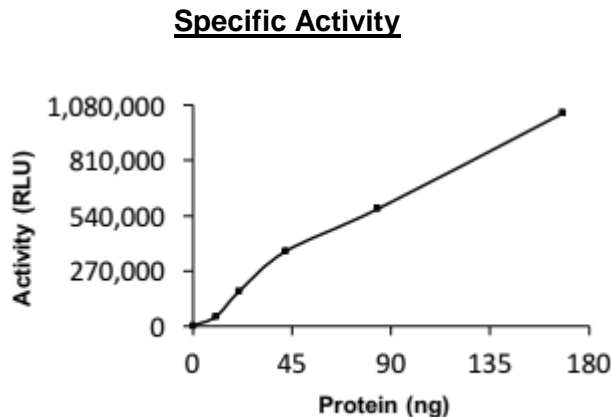
References:

1. Mondal, S., *et al.*, *Anal. Biochem.* 2016; **510**: 41-51.
2. Rape, M., *et al.*, *Nature.* 2004; **432**: 588-595.

Description: Recombinant full-length Human UBE2C, also known as dJ447F3.2, and UBCH10. Genbank Accession No. NM_007019, with an N-terminal His-tag, expressed in E. Coli. MW=21 kDa.

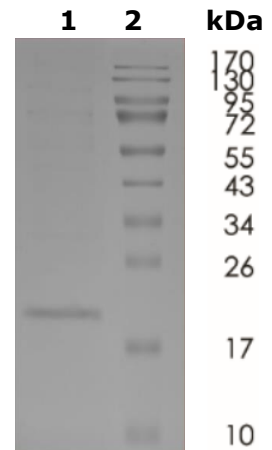
Applications: Useful for the study of ubiquitin enzyme kinetics, and as a positive control for Western blot studies.

Quality Assurance



4-20% SDS-PAGE
Coomassie staining

Lane 1:
 UBE2C
Lane 2:
 Protein Marker
MW: 21 kDa
Purity: ≥90%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com