

Data Sheet

Ubch5c, His-Tag

Human, Recombinant, N-terminal His tag
Catalog #: 80313
Lot #: 130405 **Conc.** 4.15

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 220mM Imidazole, 0.04% Tween20, 20% glycerol
Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

References:

1. Markson G. et al., Genome Res. (2009) In press.
2. Nuber U. and Scheffner M., J. Biol. Chem. 274 (11), 7576-7582 (1999).

Description: Human Ubch5c, also known as UBE2D3, (GenBank Accession No. U39318.1), amino-acids 2-end, with N-terminal His tag, MW=18 kDa, expressed in an E.Coli expression system.

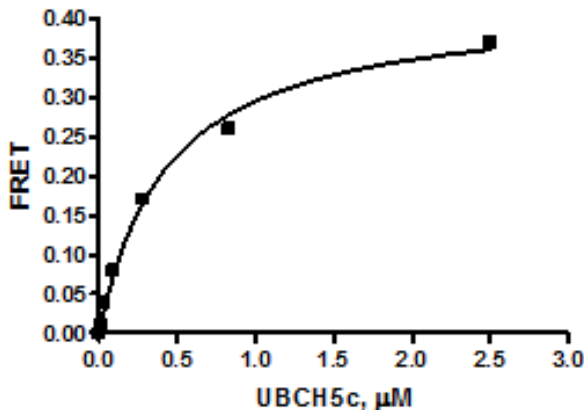
Assay Description: 10 uL of reaction mixture contains 50 mM Tris, pH 7.4, 5 mM MgCl₂, 0.01% Tween, 1 mM DTT, 0.2 μM Ub[Biotin], 50 nM Ube1, 1 mM ATP, and various amount of E2. The reaction was initiated with the addition of ATP, and incubated at 30C for 3 hours. Tb labeled anti-His antibody donor and xl665 labeled acceptor was added and incubated at room temperature for hour.

Application:

Useful in conjunction with E1 and E3 for the studies of enzyme kinetics, screening inhibitors and selectivity

Quality Assurance

Specific Activity

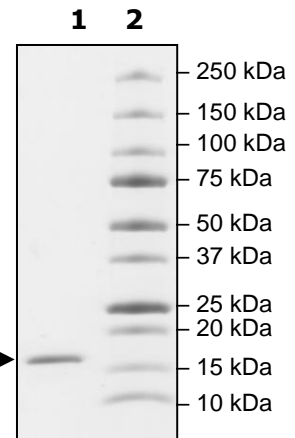


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
UBCH5C
Lane 2:
Protein Marker

Purity: $\geq 90\%$
(total)

MW: 18 kDa



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

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com