

## 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  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Protein may be diluted to  $\geq 100 \mu\text{g/ml}$  in PBS + glycerol and stored at  $-80^{\circ}\text{C}$ .

**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 (GenBank Accession No. NM\_001300795), amino-acids 2-end, with N-terminal His tag, MW=17 kDa, expressed in an E.Coli expression system.

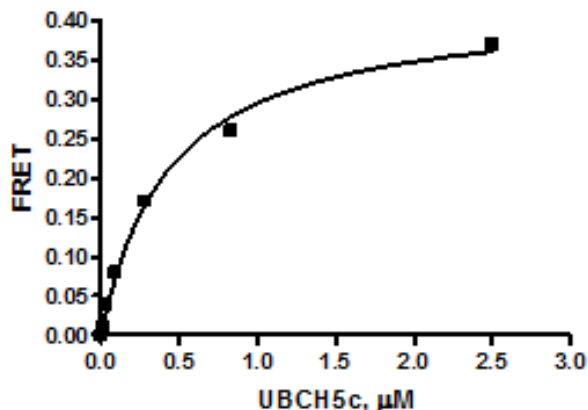
**Assay Description:** 10 uL of reaction mixture contains 50 mM Tris, pH 7.4, 5 mM MgCl<sub>2</sub>, 0.01% Tween, 1 mM DTT, 0.2  $\mu\text{M}$  Ub[Biotin], 50 nM Ube1, 1 mM ATP, and various amount of E2. The reaction was initiated with the addition of ATP, and incubated at 30C for 3 hours. Tb labeled 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 profiling.

### Quality Assurance

**Specific Activity**

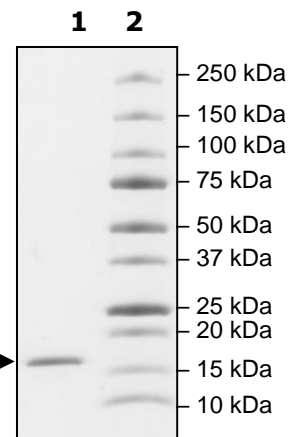


**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
UBCH5C  
**Lane 2:**  
Protein Marker

**Purity:**  $\geq 38\%$   
(total)

**MW:** 17.5 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](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)