

Data Sheet

Ubch5a, His-Tag

Human, Recombinant, N-terminal His tag

Catalog #: 80315

Lot #: 140115-B **Conc.:** 0.71 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 225 mM imidazole, 10% glycerol, and 3 mM DTT.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Protein may be diluted to ≥ 100 $\mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

References:

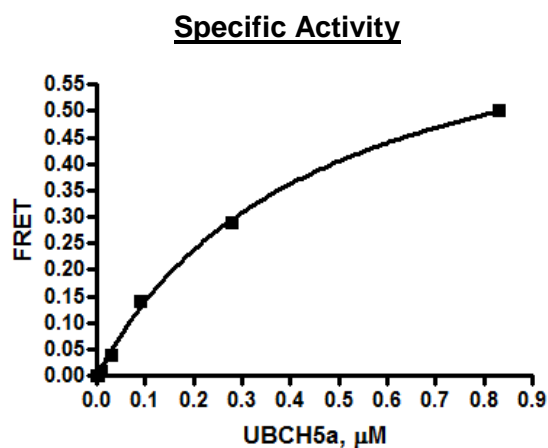
1. Markson G. et al., Genome Res. (2009) In press.
2. Nuber U. and Scheffner M., J. Biol. Chem. 274 (11), 7576-7582 (1999).

Description: Human Ubch5a (GenBank Accession No. NM_003338), amino-acids 2-end, with N-terminal His tag, MW=17 kDa, expressed in an E.Coli expression system.

Assay Conditions: 10 μL of reaction mixture contains 50 mM Tris, pH 7.4, 5 mM MgCl_2 , 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 30°C 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.

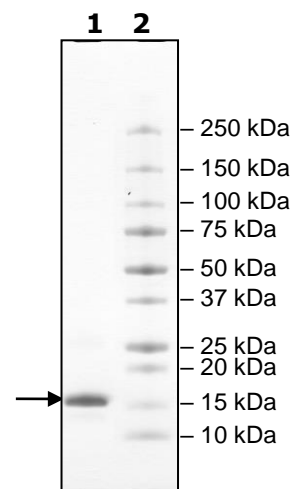
Quality Assurance



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

Lane 1:
2.70 μg
Lane 2:
Protein Marker

MW: 17 kDa
Purity: $\geq 90\%$



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

To place your order, please contact us by Phone **1.858.202.1401** 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