

Data Sheet

Ubch5b, His-Tag

Human, Recombinant, N-terminal His-tag

Catalog #: 80314

Lot#: 160707-1

Conc. 1.35 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween 20, 20% glycerol, 105 mM Imidazole.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

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).

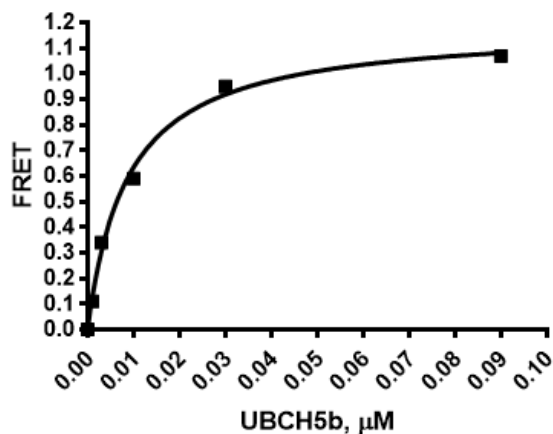
Description: Human Ubch5b also known as UBE2D2, GenBank Accession No. NM_003339, a.a 2-end, with N-terminal His-tag, expressed in an *E. coli* expression system. MW=18 kDa.

Assay Description: 10 μL of reaction mixture contains 50 mM Tris, pH 7.4, 5 mM MgCl_2 , 0.01% Tween, 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

Specific Activity



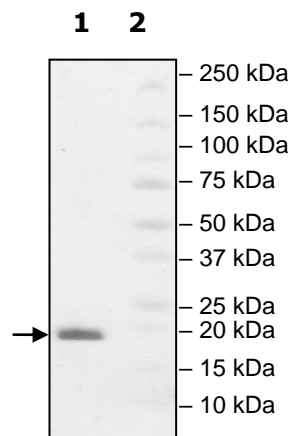
4-20% SDS-PAGE Coomassie staining

Lane 1:
UBE2D2

Lane 2:
Protein Marker

Purity: $\geq 90\%$

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, Fax 1.858.259.5734

Or you can Email us at: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com