

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).

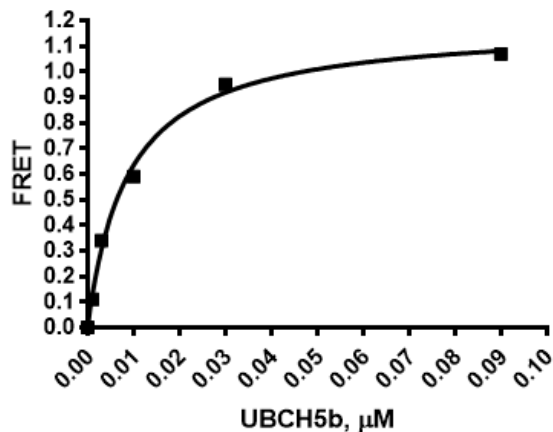
Description: Human UbcH5b also known as
 UBE2D2, GenBank Accession No.
 U39317.1, a.a 2-end, with N-terminal His-
 tag, expressed in an l expression system.
 MW=18 kDa,

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 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



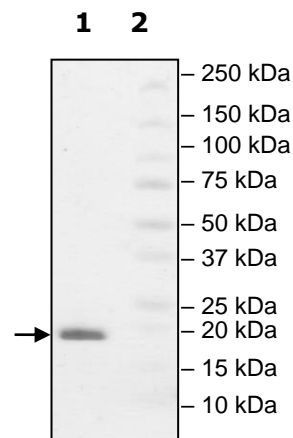
**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 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 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