

Data Sheet

Ubch13 (UBE2N)

Human, recombinant, N-terminal His-tag

Catalog #: 80323

Lot #: 91119

Conc.: 2.11 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0,
124 mM NaCl, 2.4 mM KCl, 250 mM
Imidazole, 10% glycerol and 3 mM DTT.

Stability: >6 months at -80°C

References:

1. Brusky, J., *et al.*, 2000. *Curr. Genet.* **37**: 168-174.
2. Andersen, P.L., *et al.*, 2005. *J. Cell Biol.* **170(54)** : 745-755.

Description:

Ubch13, also known as UBE2N (GenBank Accession No. NM_003348), amino acids 2-end, with N-terminal His-tag, MW = 17.6 kDa, expressed in an *E. coli* cell expression system.

Application:

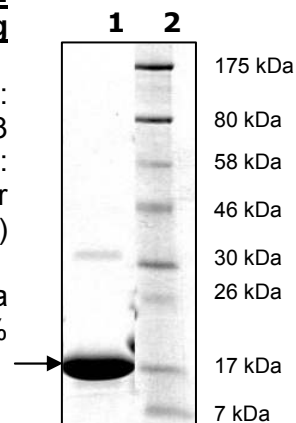
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20 % SDS-PAGE Coomassie staining

Lane 1:
42.2 µg Ubch13
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 17.6 kDa
Purity: ≥95%



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

To place your order, please contact us by Phone Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com