

Data Sheet

UBE2K (UBC1), His-Tag Protein

Human, Recombinant, with N-terminal His-tag

Catalog #: 79373

Lot #: 171023 **Conc.:** 0.1 mg/ml

Description: Recombinant full length UBE2K (UBC1) protein also known as E2-25K, HIP2, HYPG, LIG, and UB1. Genbank Accession NM_005339, with an N-terminal His-tag, expressed in E. coli cells. MW = 25kDa.

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

Formulated in: 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 25% glycerol.

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

References:

1. Kalchman, M. A. *et.al. J. Biol. Chem.* **271**: 19385-19394, 1996.
2. De Pril, R. *et al. Mol Cell Neurosci.* 2007 Jan;**34(1)**:10-9.

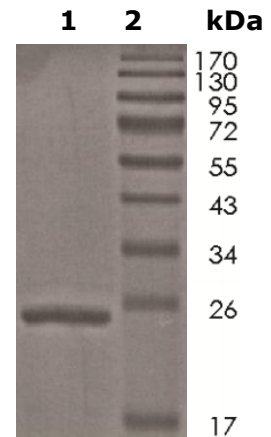
Quality Assurance

10% SDS-PAGE Coomassie staining

Lane 1:
UBE2K Protein

Lane 2:
Protein Marker

MW: ~25 kDa
Purity: ≥90%



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com