

Data Sheet

UBA6 (UBE1L2)

Human, recombinant, N-terminal Flag tag
Catalog #: 80303
Lot #: 120824 **Conc:** 0.13 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 80 ng/μl Flag
peptide, 20% Glycerol and 3 mM DTT.

Stability: >6 months at -80 °C

References:

1. Groettrup M. *et al.*, *Trends Biochem. Sci.* **33** (5), 230-237 (2008).
2. Pelzer C. *et al.*, *J. Biol. Chem.* **282** (32), 23010-23014 (2007).

Description:

Human UBA6 (UBE1L2) (GenBank Accession No. NM_018227), amino-acids 2-end, with N-terminal Flag tag, MW=118 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

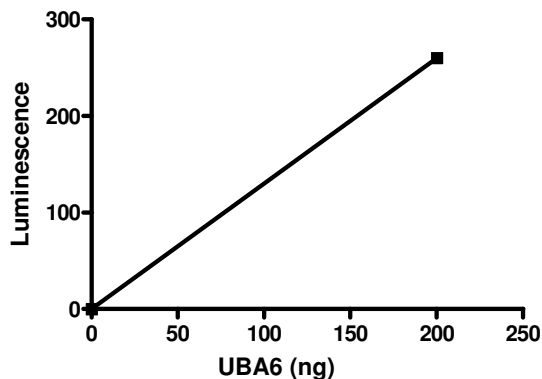
Assay Conditions: Reaction mixture containing UBA6 (UBE1L2) as a positive control, Ubc5a (500 ng), and Flag-Smurf2 (500 ng) was incubated at 37 °C for 30 minutes with biotinylated ubiquitin.

Application:

Useful in conjunction with E2 and E3 for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

Specific Activity



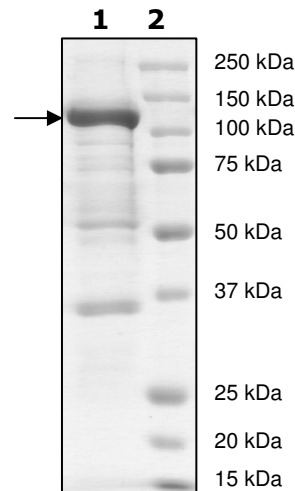
10% SDS-PAGE Coomassie staining

Lane 1:
2.6 μg UBA6

Lane 2:
Protein Marker
Biorad (#161-0373)

MW: 118 kDa

Purity: ≥57%



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