

Data Sheet

USP2

Human, recombinant, N-terminal His-tag
Catalog #: 80392
 Lot#: 120928 Conc.: 1.0 mg/ml

Formulated in: 40 mM Tris-HCl, pH 7.5,
 110 mM NaCl, 2.2 mM KCl, 15 mM β-
 mercaptoethanol, 250 mM Imidazole,
 and 20% glycerol

Stability: >6 months at -80°C

References:

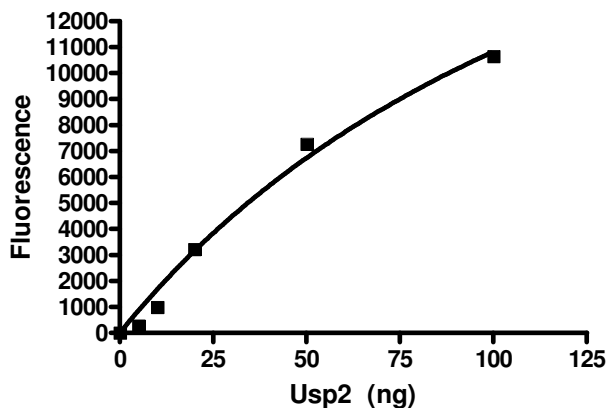
1. Stevenson, L.F., *et al.* (2010). *EMBO J.* **140(3)**: 384-96.
2. Priolo, C., *et al.* (2006). *Cancer Res.* **66(17)**: 8625-32.
3. Haimerl, F., *et al.* (2009). *J. Biol. Chem.* **284(1)**: 495-504.

Description: Human USP2 (Ubiquitin-Specific Protease 2) or UPB41, GenBank Accession No. NM_004205, amino-acids 259-605(end) with an N-terminal His-tag, MW=40.9 kDa, expressed in an *E. coli* cell expression system.

Specific Activity: 247 pmol/min/μg
 Assay Conditions: Reaction was performed in 50 mM Tris, pH 7.4, 1 mM DTT, 0.5 mM EDTA, 500 nM Ub-AMC, and USP2 in a 50 μL reaction. Reaction was incubated at 37°C for 15 min and fluorescent signal was measured at excitation = 340 nm, emission = 460 nm.

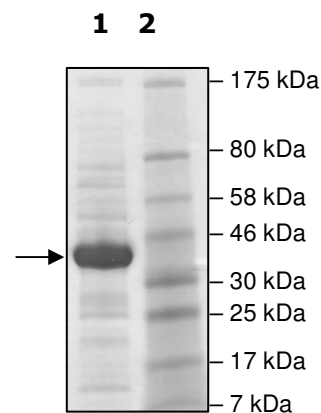
Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



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

Lane 1:
 5 μg USP2
Lane 2:
 Protein Marker
MW: 40.9 kDa
Purity: ≥85%





6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

**MHHHHHHHNSKSAQGLAGLRNLGNTCFMNSILQCLSNTRELRDYCLQRLYMRDLHHG
SNAHTALVEEFAKLIQTIWTSSPNDVVSPSEFKTQIQRYAPRFVGYNQDQAQEFRLFL
LDGLHNEVNRVTLRPKSNPENLDHLPDDEKGRQMWRKYLEREDSRIGDLFVGQLKS
SLTCTDCGYCSTVFDPFWDLSLPIAKRGYPEVTLMDCMRLFTKEDVLDGDEKPTCCR
CRGRKRCIKKFSIQRFKILVLHLKRFSESRI RTSKLTTFVNFPLRDLDLREFASENTNH
AVYNLYAVSNHSGTTMGGHYTAYCRSPGTGEWHTFNDSSVTPMSSSQVRTSDAYLL
FYELASPPSRM**

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