

Data Sheet

USP10, FLAG-tag

Human, recombinant, N-terminal FLAG-tag
Catalog #: 80360
Lot #: 140501-2 **Conc.:** 0.22 mg/ml

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

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 - 0.5%).

References:

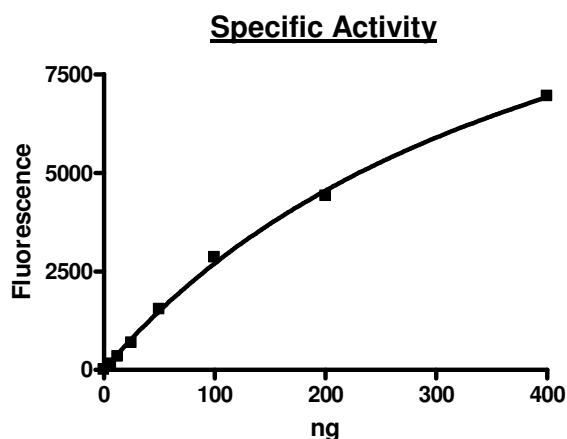
1. Yuan, J., *et al.* (2010). *Cell* **140(3)**: 384-96.
2. Soncini, C., *et al.* (2001). *Oncogene* **20(29)**:3869-79.
3. Faus, H., *et al.* (2005). *Mol. Cell. Endocrinol.* **245(1-2)**:138-46.

Description: Human USP10 (Ubiquitin-Specific Protease 10) or Ubiquitin C-terminal Hydrolase 10, GenBank Accession No. NM_005153, amino-acids 2-798(end) with an N-terminal FLAG-tag, MW=88 kDa, expressed in a baculovirus-infected Sf9 cell expression system.

Specific Activity: 1.5 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 various amount of USP10 in a volume of 50 μl. Reaction was incubated at 37°C for 30 min and fluorescent signal was measured at excitation of 360 nm, and emission at 460 nm.

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

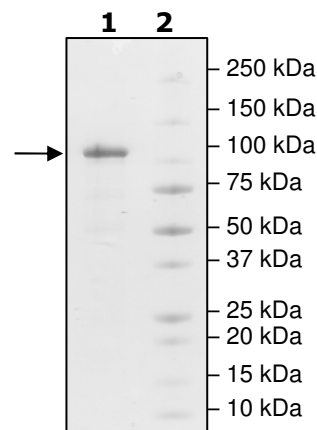
Quality Assurance



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

Lane 1:
 2.9 μg USP10
Lane 2:
 Protein Marker

MW: 88 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