

Data Sheet

USP8

Human, recombinant, N-terminal FLAG-tag

Catalog #: 80358

Lot#: 140522-A **Conc.:** 0.11 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 0.04%
 Tween20, 80 ug/ml flag peptide, 20%
 glycerol

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

References:

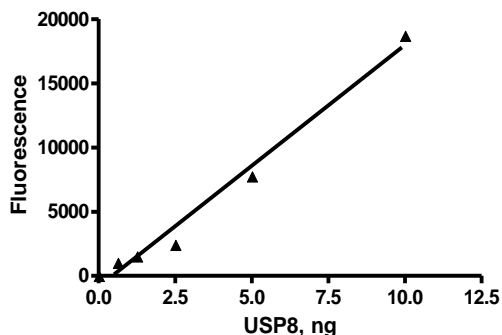
1. Naviglio, S., *et al.* (1998). *EMBO J.* **17(12)**:3241-50.
2. Niendorf, S., *et al.* (2007). *Mol Cell Biol.* **27(13)**:5029-39.
3. Mizuno, E., *et al.* (2005). *Mol. Biol. Cell.* **16(11)**: 5163-74.

Description: Human USP8, also known as Ubiquitin-Specific Protease 8, and UBPY, GenBank Accession No. NM_005154, amino-acids 2-1118(end) with an N-terminal FLAG-tag, expressed in a baculovirus-infected Sf9 cell expression system. MW=128 kDa.

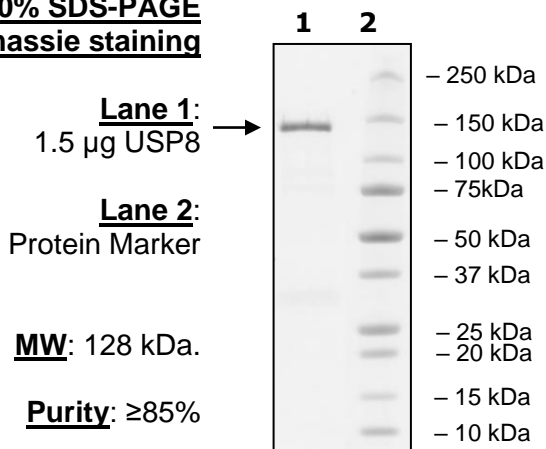
Specific Activity: 20 pmole/min/ μg
 Assay Conditions: Reaction was performed in 50 mM Tris, pH 7.4, 1 mM DTT, 0.5 mM EDTA, 500 nM Ub-AMC, and USP8 in a 50 μL reaction. Reaction was incubated at 37°C for 15 minutes and fluorescent signal was measured at excitation = 340 nm, emission = 460 nm.

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

Quality Assurance



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



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com