

Data Sheet

USP2, FLAG-Tag

Human, Recombinant, N-terminal FLAG-tag
Catalog #: 80352
Lot #: 181120-1 **Conc.:** 0.28 mg/ml

Formulated in: 40mM Tris-HCl pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 3 mM DTT,
 100ug/mL FLAG peptide 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:

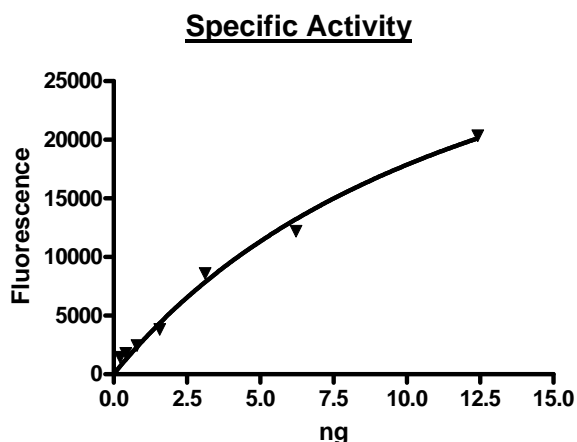
1. Stevenson, L.F., *et al.* (2010). *EMBO J.* **140(3)**: 384-96.
2. Priolo, C., *et al.* (2006). *Cancer Res.* **66(17)**: 8625-32.
3. Haimerl, F., *et al.* (2009). *J. Biol. Chem.* **284(1)**: 495-504.

Description: Human USP2 (Ubiquitin-Specific Protease 2) or UPB41, GenBank Accession No. NM_004205, amino-acids 2-605(end) with an N-terminal FLAG-tag, MW=69 kDa, expressed in a baculovirus-infected Sf9 cell expression system.

Specific Activity: 8.9 pmol/min/ μg
Assay Conditions: The enzymatic reactions were conducted in duplicate at room temperature for 30 minutes in a 50 μl mixture containing 50 mM Tris pH 7.4, 1 mM DTT, 0.5 mM EDTA, 500 nM Ub-AMC, and various amount of USP2 in a volume of 50 μl . Fluorescent signal was measured at excitation of 360 nm and emission at 460 nm after a 15 min incubation at 37°C .

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

Quality Assurance



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

Lane 1:
 2 μg USP2
Lane 2:
 Protein Marker
MW: 69 kDa
Purity: 80%

