

Data Sheet

USP5(IsoT), FLAG-tag

Human, recombinant, N-terminal FLAG-tag

Catalog #: 80355

Lot#: 120213

Conc.: 0.37 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, 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. Dayal, S., *et al.* (2009). *J. Biol. Chem.* **284(8)**:5030-41.
2. Reyes-Turcu, F.E., *et al.* (2008). *J. Biol. Chem.* **283(28)**:19581-92.

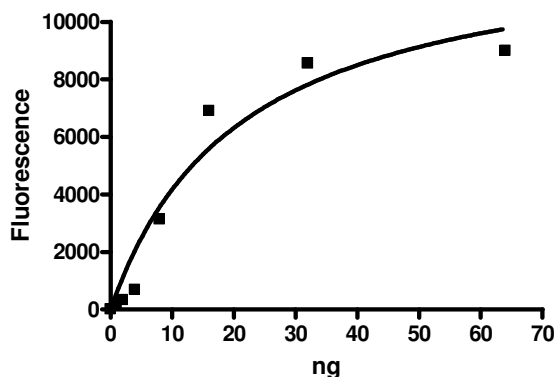
Description: Human USP5 (Ubiquitin-Specific Protease 5) or Isopeptidase T (IsoT), GenBank Accession No. NM_001098536, amino-acids 2-858(end) with an N-terminal FLAG-tag, MW=96 kDa, expressed in a baculovirus-infected Sf9 cell expression system.

Specific Activity: 12 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 USP5 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.

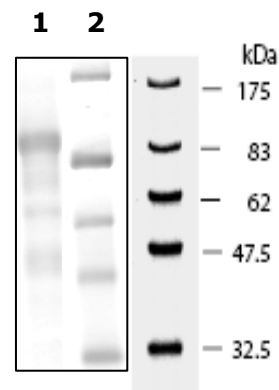
Quality Assurance



10% SDS-PAGE →
Coomassie staining

Lane 1:
2.3 μg USP5
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 96 kDa
Purity: ≥51%



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