

Data Sheet

USP14

Human, recombinant, N-terminal FLAG-tag

Catalog #: 80364

Lot#: 120227

Conc.: 0.42 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 80 ng/ul Flag
 peptide, 3mM DTT, 20% Glycerol.

Stability: >6 months at -80 °C

References:

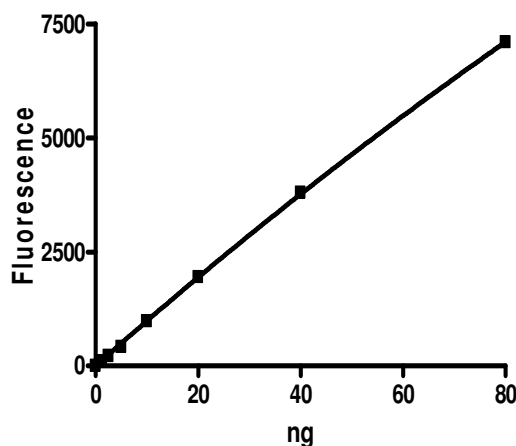
1. Chen, P.C., *et al.* (2009). *J. Neurosci.* **29(35)**:10909-19.
2. Chen, Y.C., *et al.* (2010). *RNA* **16(5)**:958-68.
3. Nagai, A., *et al.* (2009). *Biochem. Biophys. Res. Commun.* **379(4)**:995-1000.

Description: Human USP14 (Ubiquitin-Specific Protease 14) or TGT (tRNA-guanine transglycosylase), GenBank Accession No. NM_005151, amino-acids 2-494(end) with an N-terminal FLAG-tag, MW=57 kDa, expressed in a baculovirus-infected Sf9 cell expression system.

Specific Activity: 5.7 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 USP14 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.

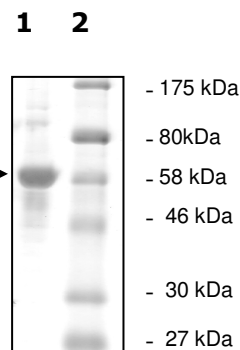
Quality Assurance



10% SDS-PAGE Coomassie staining

Lane 1:
2.8 μg USP14
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 57 kDa
Purity: ≥75%



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