

Data Sheet

UCHL3

Human, recombinant, N-terminal His-tag

Catalog #: 80353

Lot#: 100826 Conc.: 2.07 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 225 mM imidazole, 10% glycerol, and 3 mM DTT.

Stability: >6 months at -80°C

References:

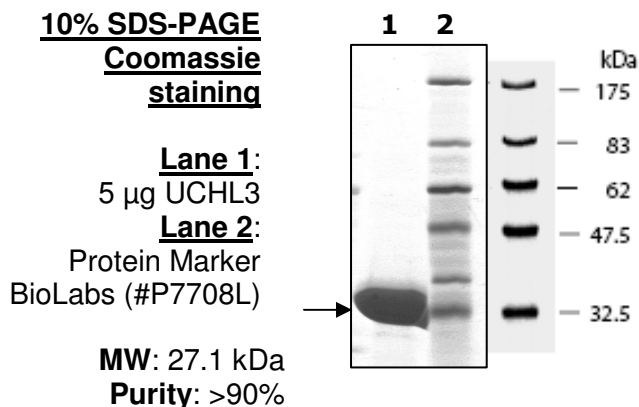
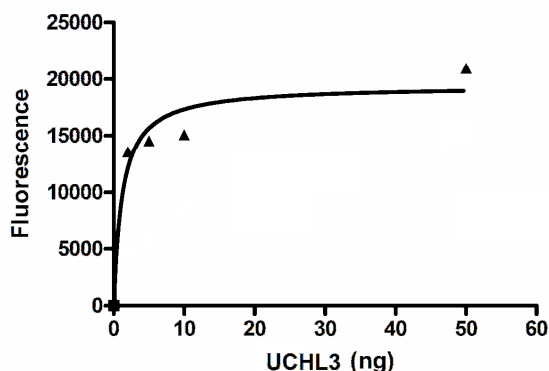
1. Wada, H., *et al.* (1998). *Biochem. Biophys. Res. Commun.* **251** (3): 688-92.
2. Nam, M.J., *et al.* (2004). *Proteomics* **3** (11): 2108-15.
3. Setsuie, R., *et al.* (2009). *Neurochem Int.* **54**(5-6):314-21.

Description: UCHL3, also known as Ubiquitin Carboxy-terminal Hydrolase L3 (GenBank Accession No. NM_006002), a.a. 2-231(end), with N-terminal His-tag, MW=27.1 kDa, expressed in *E. coli*.

Specific Activity: 20,733 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 UCHL3 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



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