

Data Sheet

UCHL1, His-Tag

Human, recombinant, N-terminal His-tag

Catalog #: 80351

Lot#: 200924-1 **Conc.:** 1.95 mg/ml

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

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

References:

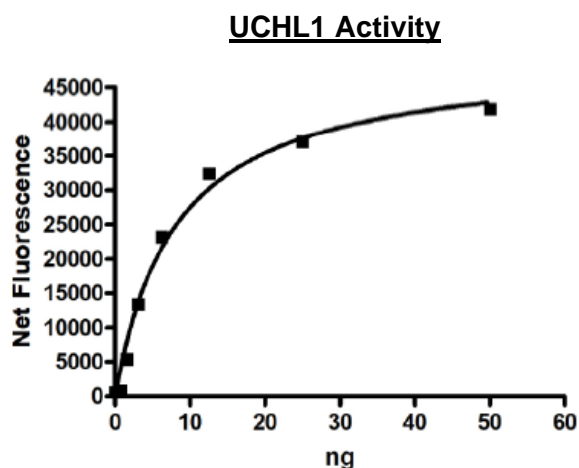
1. Gong, B., *et al.*, (2006). *Cell* **126** (4): 775–788.
2. Healy DG, *et al.*, (2005). *Cell Tissue Res.* **318** (1): 189–94.

Description: Human Ubiquitin Carboxy-terminal Hydrolase (UCHL1), also known as PGP9.5, GenBank Accession No. NM_004181, a.a. 2-223, with N-terminal His-tag, expressed in *E. coli*. expression system. MW=26 kDa.

Specific Activity: 6000 pmol/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 UCHL1. Reaction was incubated at 37°C for 15 min. and fluorescent signal was measured at excitation of 340 nm, and emission at 460 nm.

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

Quality Assurance

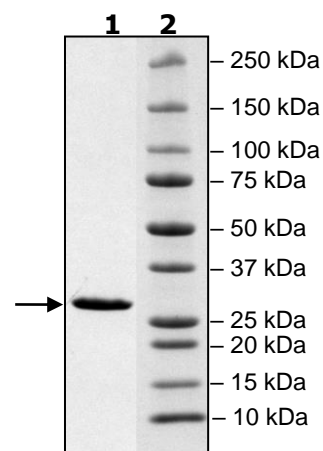


**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 2 μg UCHL1
Lane 2:
 Protein Marker

MW: 26 kDa

Purity: >90%



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