

UCHL3 (Ubiquitin C-terminal hydrolase L3)

Cat. # DB101

Background: UCHL3 have 52% amino acid identity with UCHL1 and is uniformly expressed in all tissues, including brain¹. The activity of UCHL3 is more than 200-fold higher than UCH-L1 when a fluorogenic ubiquitin substrate is used².

Alternate names: Ubiquitin carboxyl-terminal hydrolase isozyme L3, Ubiquitin thioesterase L3, UCH-L3

Product Information

Molecular Weight:	25kDa
Quantity:	25µg
Physical State:	Liquid
Source:	Human Recombinant
Activity:	This enzyme cleaves Ub-AMC
Storage:	-80°C. Avoid repeated freeze/thaw cycles

References

1. Mayer, A. N. and Wilkinson, K. D. *Detection, resolution, and nomenclature of multiple ubiquitin carboxyl-terminal esterases from bovine calf thymus*. *Biochemistry* 1989. 28: p166-172
2. Larsen C. et al. *Substrate specificity of deubiquitinating enzymes: ubiquitin C-terminal hydrolases*. *Biochemistry* 1998. 37: p3358–3368.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved