

UCH-L3 (Ubiquitin C-terminal hydrolase L3)

Cat. # DB101

Background

UCH-L3 is a deubiquitinase that controls levels of cellular ubiquitin through the processing of ubiquitin precursors and ubiquitinated proteins. It is a thiol protease that recognizes and hydrolyzes a peptide bond at the C-terminal glycine of either ubiquitin or NEDD8. The activity of UCH-L3 is more than 200-fold higher than that of UCH-L1 when a fluorogenic form of ubiquitin is used as a substrate.

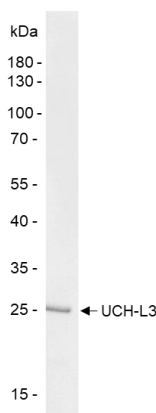
Alternate names

Ubiquitin carboxyl-terminal hydrolase isozyme L3, Ubiquitin thioesterase L3, UCH-L3

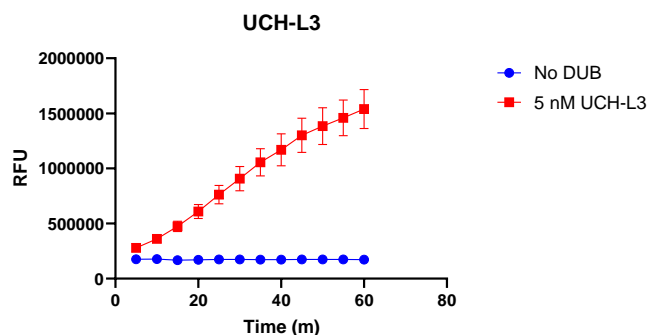
Product Information

Purity	≥ 95% by SDS-PAGE
Molecular Weight	25 kDa
Quantity	25 µg
Physical State	Liquid
Source	Human Recombinant
Tag	None
Activity	This enzyme cleaves Ub-AMC
Storage	-80°C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified UCH-L3. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of UCH-L3. 5 nM UCH-L3 was tested in a Ub-Rh110 assay showing robust DUB activity.

References

- Hafez, N., et al., Eur. J Med Chem., 2022. 227:113970.
- Lee. J E., et al., Anim Cells Sys (Seoul), 2021. 25(5):312-322.

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