

UCH-L5 (Ubiquitin C-terminal hydrolase L5)

Cat. # DB102

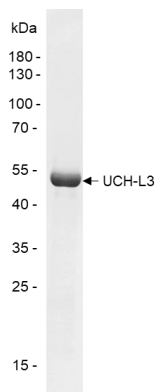
Background UCH-L5, or Ubiquitin C-Terminal Hydrolase L5, is a deubiquitinating enzyme that plays a vital role in the regulation of protein degradation within cells. As a member of the ubiquitin carboxyl-terminal hydrolase (UCH) family, UCH-L5 functions to remove ubiquitin molecules from protein substrates, thereby preventing their degradation via the proteasome or lysosome. UCH-L5 is particularly known for its role in the 19S proteasome, where it acts as a deubiquitinating enzyme to process and recycle ubiquitin from substrate proteins before their proteolytic degradation. This activity is crucial for maintaining protein quality control and ensuring proper protein turnover within cells.

Alternate names AD-019, CGI-70, Ubiquitin carboxyl-terminal hydrolase isozyme L5, Ubiquitin C-terminal hydrolase UCH37, Ubiquitin thioesterase L5, UCH37, UCH-L5

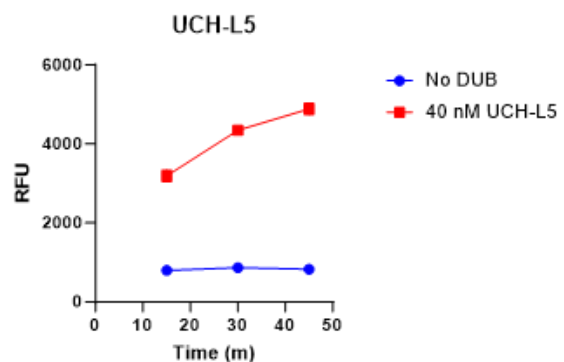
Product Information

Purity	≥ 95% by SDS-PAGE
Molecular Weight	49 kDa
Quantity	25 µg
Physical State	Liquid, 50mM Tris, pH8.0, 150mM NaCl, 10% glycerol
Source	Human Recombinant
Tag	His6-SUMO
Activity	This enzyme is active in the Ub-CHOP assay or with ubiquitin-AML (Cat. # SI260)
Storage	-80° C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified UCH-L5. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of UCH-L5. 40 nM UCH-L5 was tested in a Ub-Rh110 assay showing robust DUB activity.

References

1. Ge, J., et al., Oncotarget., 2017. 8(69):113635-113649.
2. Sharma, A., et al., Int J Mol Sci., 2020. 21(11):3910.

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