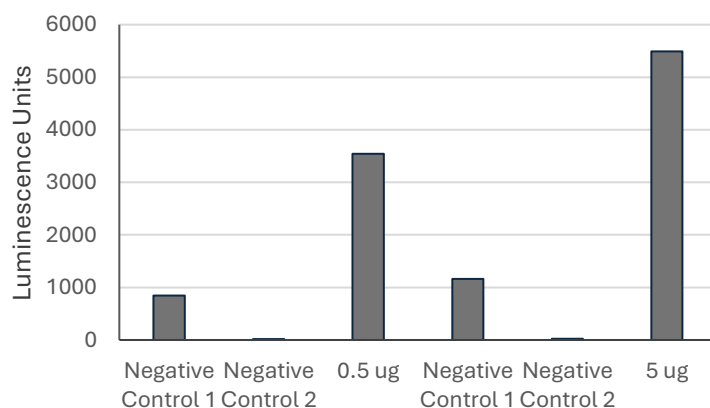


Background	UCHL5 is a member of small family of DUBs referred to as the Ub C-terminal hydrolases (UCHs). UCHL5 is the only UCH family protease, that is associated with mammalian proteasomes. It is responsible for the ubiquitin isopeptidase activity in the PA700 (19S) proteasome regulatory complex. UCHL5 exclusively cleaves branched Ub chains. K48 linkages are readily removed in branched chains.
Alternate Names	AD-019, CGI-70, Ubiquitin carboxyl-terminal hydrolase isozyme L5, Ubiquitin C-terminal hydrolase UCH37, Ubiquitin thioesterase L5, UCH37, UCH-L5
Application(s)	ELISA, Western Blot (Recombinant protein was detected at a final concentration of 10.14µg/mL).

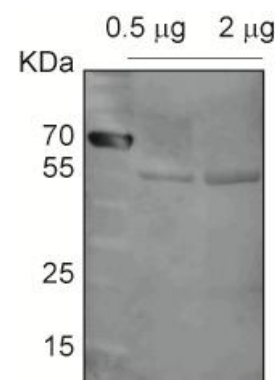
Product Specifications

Purity	≥ 90% estimated by SDS-PAGE
Molecular Weight	49 kDa
Quantity	250 µg
Physical State	Liquid. Supplied in phosphate buffered saline pH 7.4.
Source & Clonality	Chicken. Polyclonal antibody to UCHL5
Species Reactivity	Human, Chicken
Immunogen	Full-length human UCHL5
Isotype	IgY
Stability & Storage	Over 1 year at -20° C. Avoid repeated freeze/thaw cycles

Product QC



Analysis of UCH-L5 using ELISA plate. UCH-L5 of 0.5 or 5ug was detected by 2.03 µg/mL of AB102 using traditional ELISA detection reagent. 2° Antibody: α-Chicken HRP (1:5000). Negative Control 1: No AB102; Negative Control 2: no detection reagent 2: no detection reagent. AB102.



SDS-Page Analysis of purified UCH-L5. Indicated amount loaded on SDS-PAGE gel followed by Western Blot using 10.14 µg/mL of UCH-L5. 2° Antibody: α-Chicken HRP (1:5000).

References

1. Deol KK., et al., Mol Cell. 2020 ;80(5) :796-809.e9.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved