

CYLD (human) polyclonal antibody

The tumor suppressor CYLD is a deubiquitylating enzyme (DUB) that antagonises NF- κ B and JNK signaling by disassembly of Lys⁶³-linked ubiquitin chains synthesised in response to cytokine stimulation. The DUB activity of CYLD is mediated by its C-terminal ubiquitin specific protease (USP) domain, pathogenic truncations and mutations of which are associated with the hypertrophic skin tumor cylindromatosis through disruption of the USP domain and subsequent loss of CYLD catalytic activity.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-PW0760-0025	25 μ l
-----------------	------------

Manuals, SDS & CofA

[View Online »](#)

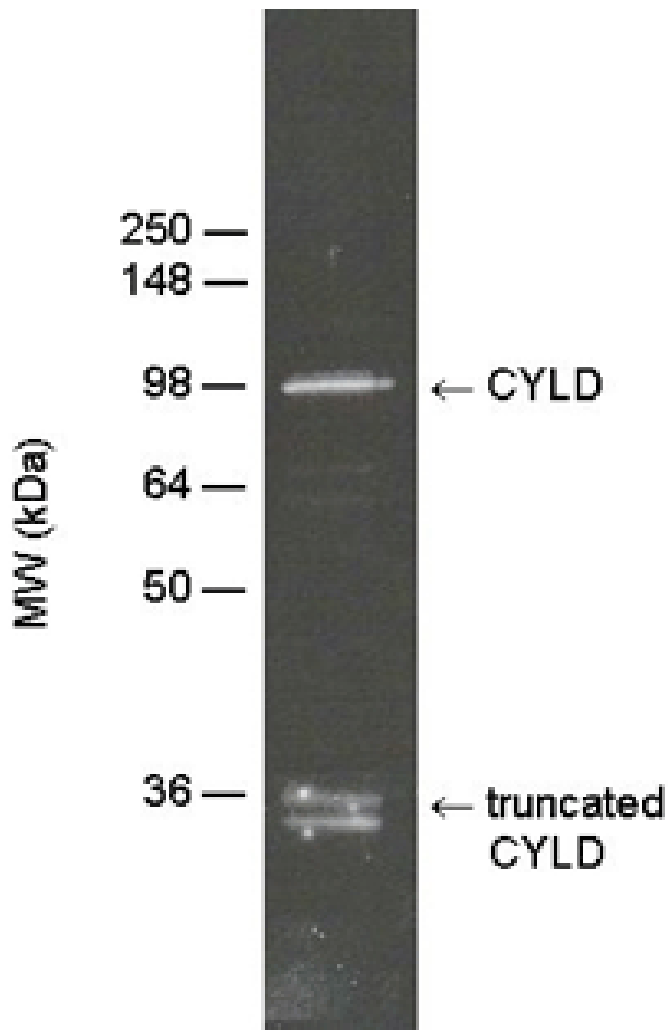


Figure: Detection of CYLD in HeLa S100 cytosolic fraction (Prod. No. BML-SW8750, 10 μ g) by Western blotting using PAb to CYLD (human) (Prod. No. BML-PW0760) at 1:5000 dilution.

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C. Dilute with PBS, pH 7.2-7.4, containing 1% normal goat serum (if a goat anti-rabbit IgG linker antibody is to be used). Store diluted antibody at +4°C, do not freeze and use within 1 month.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Probable ubiquitin carboxyl-terminal hydrolase CYLD
Application	WB
Formulation	Liquid. In PBS containing 0.01% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 336-353 of human CYLD.
Purity Detail	Purified by salt precipitation.
Recommendation Dilutions/Conditions	Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Human
UniProt ID	Q9NQC7
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)