

# Huntingtin polyclonal antibody

Huntington's disease (HD) is a neurodegenerative disorder caused by an expanding polyglutamine repeat in the huntingtin gene (>35 repeats results in HD onset). Expansion of this polyglutamine repeat may induce a toxic gain of function perhaps through interactions with other cellular proteins. The huntingtin protein is thought to be necessary for neuronal survival and is involved in synaptic vesicle trafficking, microtubule binding and may also have a role in apoptosis. SUMO modification at the N-terminus of the huntingtin protein is also believed to play a role in the neurodegenerative process.

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

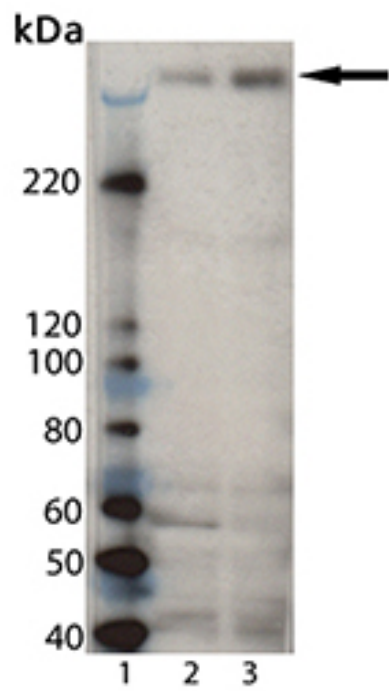
## Ordering Information

[Order Online »](#)

BML-PW0595A-0025	25µg
BML-PW0595A-0100	100µg

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Huntingtin pAb (Prod. No. BML-PW0595A): Lane 1: MW Marker, Lane 2: Mouse Brain Tissue Extract (Prod. No. ADI-LYT-MB100), Lane 3: Rat Brain Tissue Extract (Prod. No. ADI-LYT-RB100).

# Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Htt, HD Protein
Application	WB
Application Notes	Detects a band of ~350kDa by Western blot.
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 2-17 of human huntingtin.
Purity Detail	Protein A affinity purified.
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	Mouse, Rat
UniProt ID	P42858
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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)