

HDAC5 polyclonal antibody

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

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

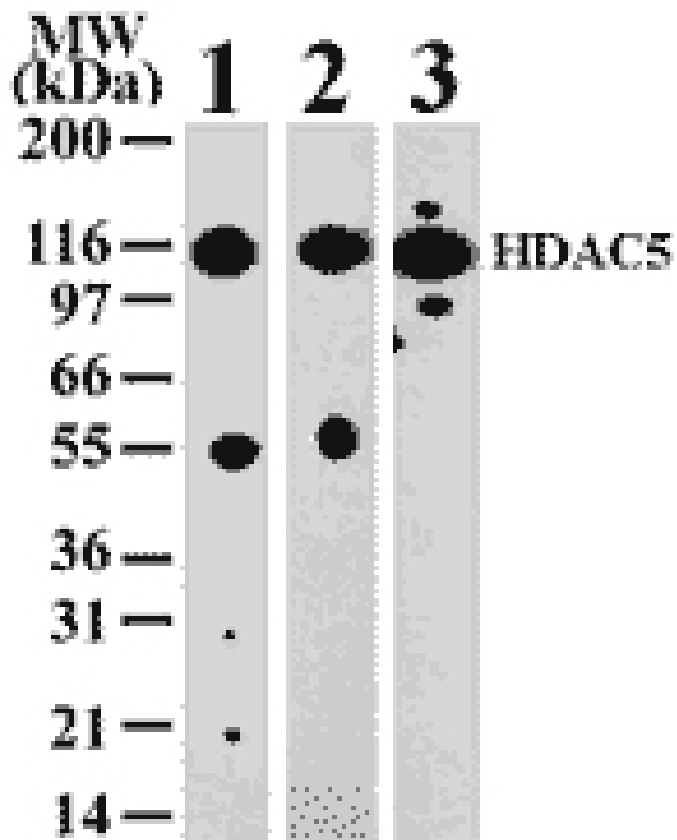


Figure: Western blot analysis of HDAC5 in 293 (lane 1), Jurkat (lane 2) and NIH-3T3 (lane 3) cell lysate using PAb to HDAC5 (Prod. No. [ALX-210-340](#)).^{*****},HDAC5 polyclonal antibody Western blot,HDAC5 polyclonal antibody Western blot,yes

ALX-210-341,210-341.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-341.gif>,**Figure:** Western blot analysis of HDAC6 in NIH-3T3 cell lysate using PAb to Histone Deacetylase 6 (Prod. No. [ALX-210-341](#)).^{*****}

Handling & Storage

Use/Stability	Stable for 6 months when stored at +4°C.
Handling	Avoid freeze/thaw cycles. After reconstitution, 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	Histone deacetylase 5
Application	ChIP, IP, WB
Application Notes	Detects a band of ~120kDa.
Formulation	Liquid. In PBS containing 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 572-587 of human HDAC5 (histone deacetylase 5).
Positive Control	NIH 3T3, 293 or Jurkat cells are recommend as positive control.
Purity Detail	Protein G-affinity purified IgG.
Recommendation Dilutions/Conditions	Western blot (2µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
UniProt ID	Q9UQL6

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)