HDAC4 (NT) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ALX-210-339-C100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

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 4

Application WB

Application NotesDetects a band of ~140kDa by Western blot.

Formulation Liquid. Purified IgG in PBS containing 0.05% BSA

and 0.05% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 194-209 of human

HDAC4 (histone deacetylase 4).

Recommendation Dilutions/Conditions Western Blot (0.5-2µg/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 P56524

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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com