

Perforin monoclonal antibody (CE2.10)

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

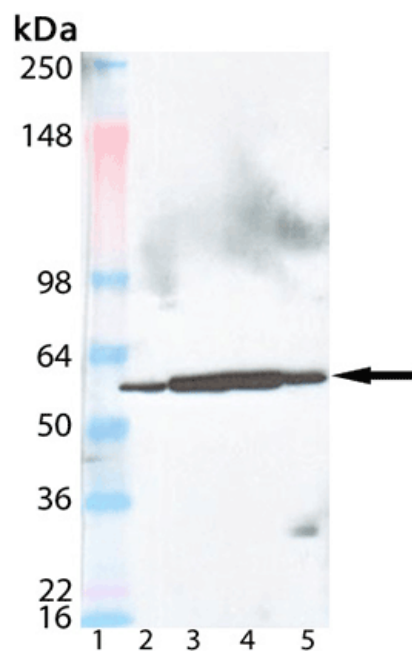
Ordering Information

[Order Online »](#)

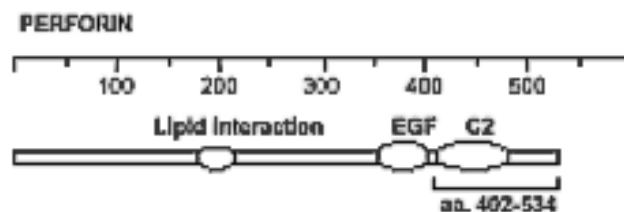
ALX-804-135-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of ALX-804-135: Lane 1: MW Marker, Lane 2: HeLa lysate, Lane 3: Jurkat lysate, Lane 4: MOLT-4 lysate, Lane 5: rat thymus lysate.



Schematic structure of mouse perforin (recombinant, aa 98-534). The antibody recognizes an epitope within aa 402-534.

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at +4°C undiluted.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	IP, WB
Application Notes	Immunoprecipitation: human and mouse Western Blot: mouse only
Clone	CE2.10
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant mouse perforin (aa 98-534).
Isotype	IgG2a
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human, Mouse
Specificity	Recognizes an epitope in the region aa 402-534.
UniProt ID	P10820
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-us@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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