

Perforin (mouse) monoclonal antibody (CB5.4)

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

Citations: 21

[View Online »](#)

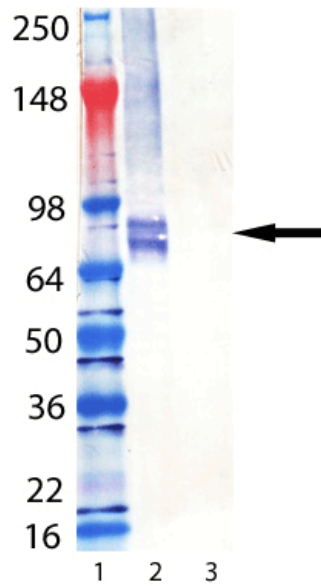
Ordering Information

[Order Online »](#)

ALX-804-057-C100	Purified - 100µg
------------------	------------------

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of Perforin (mouse) mAb (CB5.4) (Prod. No. ALX-804-057): Lane 1: MW Marker, Lane 2: CTLL-2 cell lysate, Lane 3: 3T3 cell lysate. Western Blot analysis of Perforin 1 expression in CTLL-2 cell lysate and absence of perforin in the mouse fibroblast cell line 3T3.

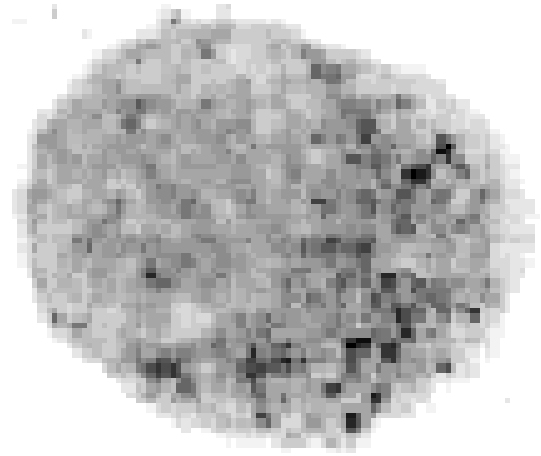
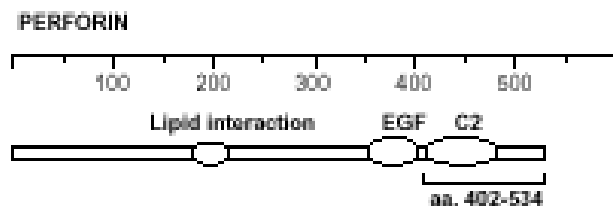


Figure: Detection of perforin in an individual mouse T cell. Perforin is located in cytoplasmic granules.

Method: A mouse T cell clone was plated onto polylysine treated glass slides, fixed and permeabilized in methanol at -20°C for 5 min, then in acetone at -20°C for 30 sec. After 3 washings in PBS, 0.1% BSA, slides were incubated with 5 $\mu\text{g}/\text{ml}$ of CB5.4 antibody, washed in PBS, incubated in anti-rat IgG-HRP, washed in PBS, and developed in appropriate substrate solution (DAB or AEC).



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

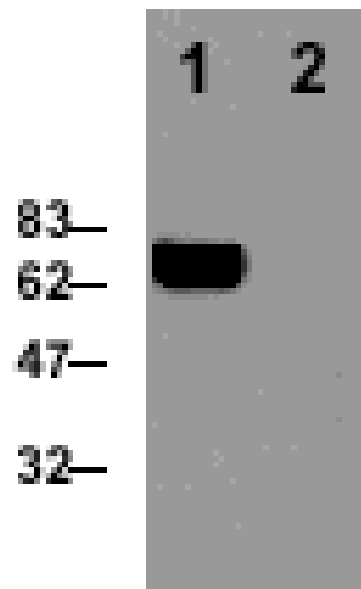


Figure: Detection of perforin in the mouse T cell clone B6.1. (lane 1). Absence of perforin in the mouse fibroblast cell line NIH.3T3 (lane 2). The protein migrates as a 66 kDa species.

Method: Cell extracts from the T cell clone B6.1 (2×10^6) were resolved by SDS-PAGE under reducing conditions, transferred to nitrocellulose and probed with the CB5.4 antibody at 1 $\mu\text{g/ml}$. Proteins were visualized using a peroxidase-conjugated antibody to rat IgG and a chemiluminescence detection system.

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at +4°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	ICC, IHC (FS), IHC (PS), IP, WB
Clone	CB5.4
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant mouse perforin (aa 98-534).
Isotype	IgG2a
Source	Purified from ascites.
Species Reactivity	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 .



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)