

# Perforin (mouse) monoclonal antibody (CB5.4)

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

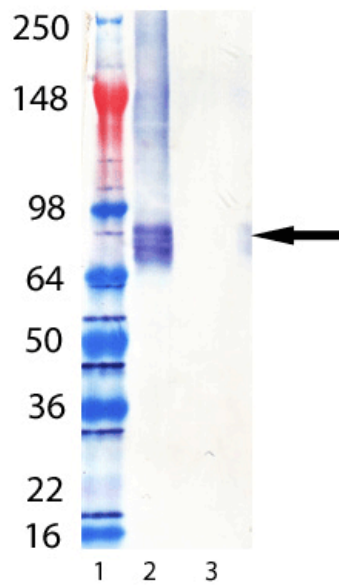
## Ordering Information

[Order Online »](#)

ENZ-ABS708-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of Perforin (mouse) mAb (CB5.4) (Prod. No. ENZ-ABS708): 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.

## Handling & Storage

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
Application Notes	Detects bands of ~66kDa and higher (depending on PTM) by Western blot.
Clone	CB5.4
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant mouse perforin (aa 98-534).
Isotype	IgG2a
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Mouse
Specificity	Recognizes an epitope in the region aa 402-534.

UniProt ID

P10820



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  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-eu@enzolifesciences.com](mailto:info-eu@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)

This antibody is covered by our **Worry-Free Guarantee**.