

# Granzyme B (human) monoclonal antibody (B18.1)

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

Citations: 6

[View Online »](#)

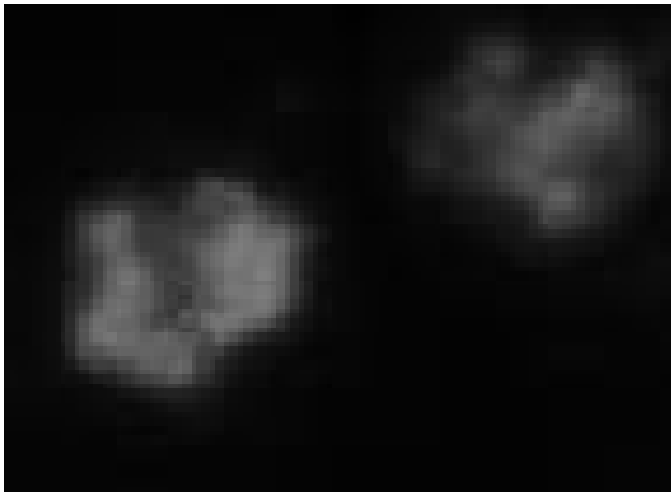
## Ordering Information

[Order Online »](#)

ALX-804-121-C100	Purified - 100µg
------------------	------------------

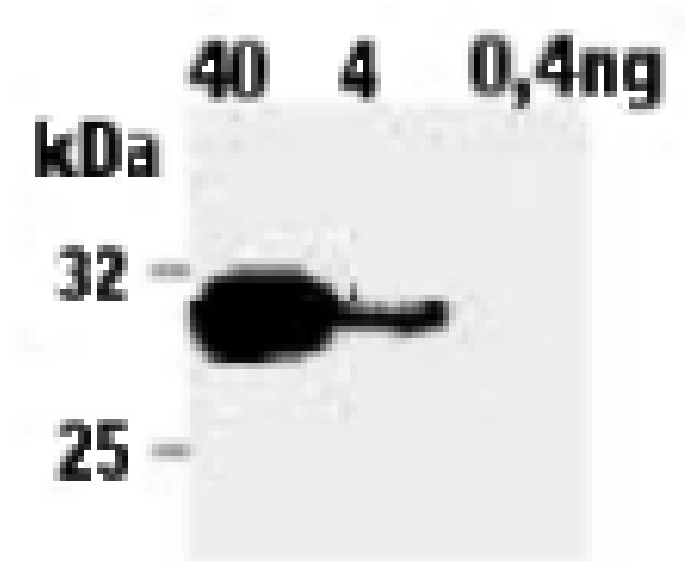
Manuals, SDS & CofA

[View Online »](#)



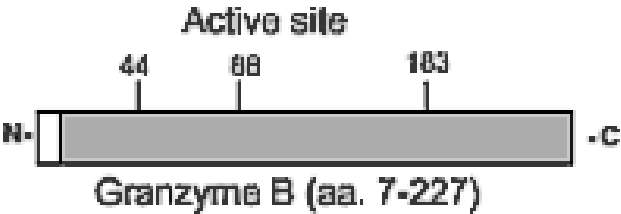
**Figure:** Detection of granzyme B in a human CD8+ T cell clone.

**Method:** Human T cells were 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 washes in PBS, 0.1% BSA, slides were incubated for 1 hr at room temperature (RT) with 20 µg/ml of B18.4 antibody in PBS, 0.1% BSA. After rinsing in PBS, a FITC-conjugated anti-mouse IgG was added for 30 min., slides washed again in PBS and visualized using a fluorescence microscope.



**Figure:** Detection of recombinant granzyme B. Granzyme B migrates as a 27 kDa species.

**Method:** Recombinant granzyme B was resolved by SDS-PAGE under reducing conditions, transferred to nitrocellulose and probed with the monoclonal anti-granzyme B antibody B18.1 at 1µg/ml. Proteins were visualized using a peroxidase-conjugated antibody to mouse IgG (SBA) and a chemiluminescence detection system.



Schematic structure of human granzyme B (recombinant, aa. 7-227).

## 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	-20°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application	Flow Cytometry, ICC, IHC (FS), IHC (PS), IP, WB
Clone	B18.1
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human granzyme B (aa 7-227).
Isotype	IgG1
Source	Purified from mouse ascites.
Species Reactivity	Human
UniProt ID	P10144
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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-  
fr@enzolifesciences.com](mailto:info-fr@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)