

# Toll-like receptor 5 monoclonal antibody (85B152.5) (R-PE conjugate)

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

Citations: 4

[View Online »](#)

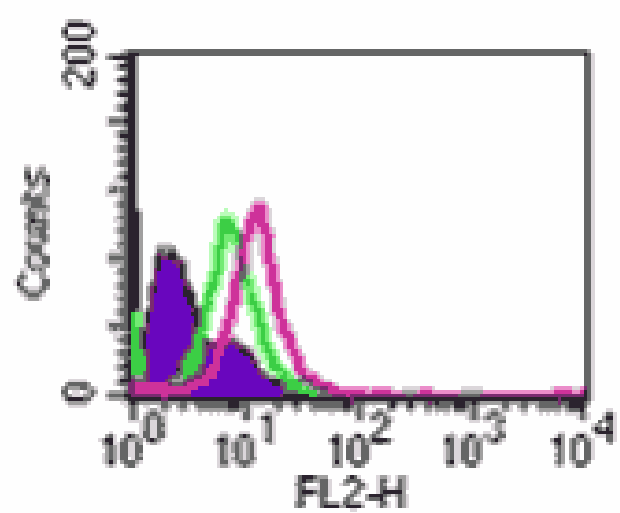
## Ordering Information

[Order Online »](#)

ALX-804-597R-C100	R-PE - 100µg
-------------------	--------------

Manuals, SDS & CofA

[View Online »](#)



# Handling & Storage

Use/Stability	Stable for 6 months when stored at +4°C.
Handling	Do not freeze. Protect from light.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	TLR5, Toll/Interleukin-1 receptor-like gene 3, TIL3
Application	Flow Cytometry
Application Notes	Flow Cytometry: Suitable for cell surface and intracellular FC.
Clone	85B152.5
Formulation	Liquid. In PBS containing 0.05% BSA and 0.05% sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to a portion of human TLR5 (Toll-like receptor 5).
Isotype	IgG2a
Positive Control	Ramos cells can be used as positive control.
Purity Detail	Protein G-affinity purified.
Species Reactivity	Dog, Human, Mouse
UniProt ID	O60602

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-  
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)