

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

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

Citations: 4

[View Online »](#)

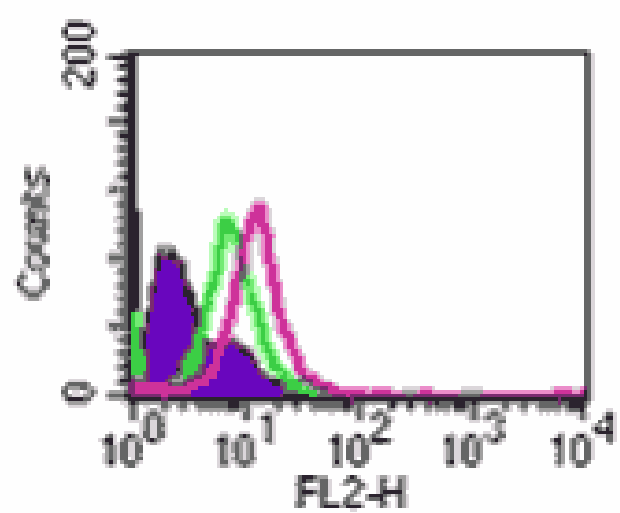
Ordering Information

[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



ENZO LIFE SCIENCES,
INC.
Phone: +1 940.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
ENZO LIFE SCIENCES
Phone: +32 3 466 0426
info-be@enzolifesciences.com

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

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

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

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