

# Toll-like receptor 5 (human) polyclonal antibody

Note: Human Toll-like Receptor 5 (TLR5) is also described as Toll/Interleukin-1 Receptor-like Gene 3 (TIL3).

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

## Ordering Information

[Order Online »](#)

ALX-210-853-C200	200µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	For maximum product recovery after thawing, centrifuge the vial before opening the cap. Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	TLR5, Toll/Interleukin-1 receptor-like gene 3, TIL3
Application	ELISA, Flow Cytometry, IHC, WB
Formulation	Affinity purified antibody in 10 mM KHPO <sub>4</sub> , 140 mM NaCl, BSA @ 1mg/ml and 0.1% sodium azide.
Host	Goat
Immunogen	Synthetic peptide corresponding to aa 151-181 (D151LSKNQIRSLYLHPSFGKLSIDFSSNQ181) of human TLR5 (Toll-like receptor 5).
Purity Detail	Epitope-affinity purified IgG.
Quality Control	Immunocytochemistry on human peripheral blood leukocytes.
Recommendation Dilutions/Conditions	ELISA (1:37,500)Western Blotting (1:50)IHC (1:250)Flow Cytometry (1:200)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	O60602

### Worry-free Guarantee



ENZO LIFE SCIENCES  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

ENZO LIFE SCIENCES  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

ENZO LIFE SCIENCES  
France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

ENZO LIFE SCIENCES  
Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

ENZO LIFE SCIENCES  
UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)