

# Toll-like receptor 10 (human) monoclonal antibody (3C10C5) (R- PE conjugate)

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

## Ordering Information

[Order Online »](#)

ALX-804-610R-T100	100 tests
-------------------	-----------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	Stable for at least 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	TLR10, CD290
Application	Flow Cytometry
Clone	3C10C5
Formulation	Liquid. In PBS containing 0.05% BSA, 0.05% sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to aa 20-474 of the extracellular domain of human TLR10.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation	Flow Cytometry (cell surface, 10 µl/106 cells (1 µg))
Dilutions/Conditions	
Species Reactivity	Human
UniProt ID	Q9BXR5

### 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  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

ENZO LIFE SCIENCES  
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)