

# Toll-like receptor 8 monoclonal antibody (44C143) (R-PE conjugate)

The Toll-like receptor (TLR) family in mammal comprises a family of transmembrane proteins characterized by multiple copies of leucine rich repeats in the extracellular domain and IL-1 receptor motif in the cytoplasmic domain. Like its counterparts in *Drosophila*, TLRs signal through adaptor molecules. The TRL family is a phylogenetically conserved mediator of innate immunity that is essential for microbial recognition. Ten human homologs of TLRs (TLR1-10) have been described. TLR8 gene contains three exons, two of which have coding function. TLR8 cDNA codes for a protein of approximate molecular weight of 120kDa.

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

Citations: 2

[View Online »](#)

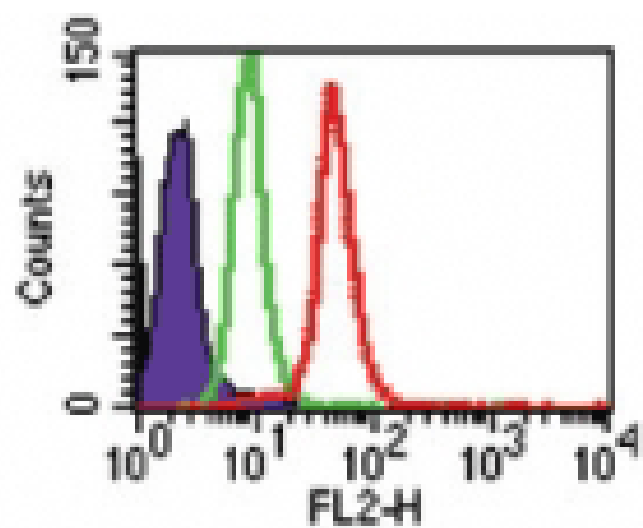
## Ordering Information

[Order Online »](#)

ALX-804-376R-C100	R-PE - 100µg
-------------------	--------------

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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	TLR8, CD288
Application	Flow Cytometry
Application Notes	Detects a band of ~120kDa by Western blot.
Clone	44C143
Formulation	Liquid. In PBS containing 0.05% BSA and 0.05% sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to aa 750-850 of the cytoplasmic portion of human TLR8 (Toll-like receptor 8).
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (intracellular, 0.5-2µg per 10 <sup>6</sup> cells; cell surface: see Xu et al. (2006) for details)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
Technical Info / Product Notes	Replacement for <b>ADI-CSA-825</b>
UniProt ID	Q9NR97

Worry-free Guarantee

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



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)