

# 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.

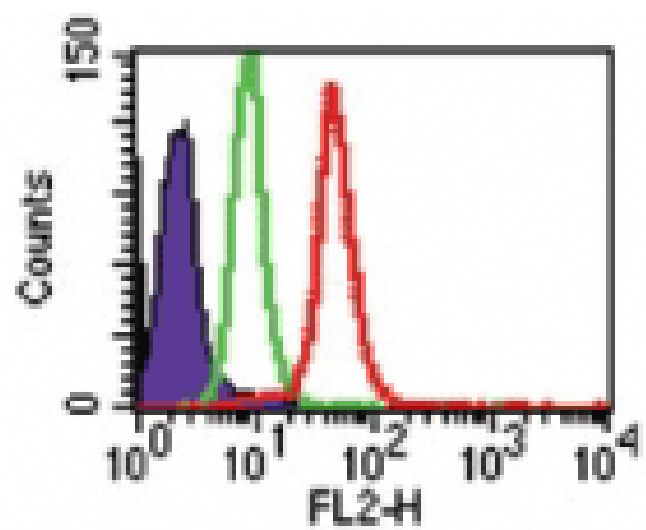
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

[View Online »](#)

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](#)

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)