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

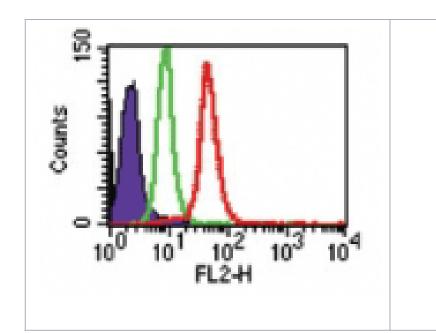
Order Online »

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 NotesDetects 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).

lsotype lgG1

Purity Detail Protein G-affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (intracellular, 0.5-2µg per 106 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 ADI-CSA-825

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com