Toll-like receptor 6 (human) monoclonal antibody (86B1153.2) (FITC conjugate)

This antibody is covered by our Worry-Free Guarantee.

Citations: 4

View Online »

Ordering Information

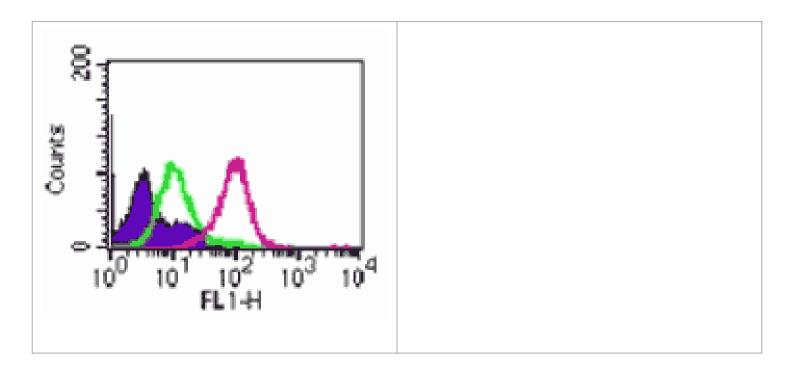
Order Online »

ALX-804-363F-C100

FITC - 100µg

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 TLR6, CD286

Application Flow Cytometry

Application Notes Flow Cytometry: Suitable for cell surface and intracellular

FC.

Clone 86B1153.2

Formulation Liquid. In 10mM TRIS containing 150mM sodium chloride

and 0.05% sodium azide.

Host Mouse

Immunogen Synthetic peptide corresponding to aa 408-424 in the

cytoplasmic portion of human TLR6 (Toll-like receptor 6).

lgG1

Positive Control Ramos cells can be used as positive control.

Purity Detail Protein G-affinity purified.

Species Reactivity Human

UniProt ID Q9Y2C9

Last modified: May 29, 2024



info-

eu@enzolifesciences.com