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.

lsotype lgG1

Purity Detail Protein G-affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (cell surface, 10 μl/106 cells (1 μg))

Species Reactivity Human

UniProt ID Q9BXR5

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com