IL-4 Recombinant monoclonal antibody (11B11) (Rat IgG1κ)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS549-0200

200µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.

Handling Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-4, B-cell IgG differentiation factor, B-cell growth

factor 1, B-cell stimulatory factor 1, BSF-1

Application ELISA, Flow Cytometry, IHC, in vivo Assay

Clone 11B11

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Rat

Immunogen BSF-1 / IL-4 purified from serum-free supernatants of

cultured EL-4 cells (a C57BL/6 thymoma) induced with 4β-

phorbol 12β-myristate 12α-acetate for 48h.

lgG1κ

Purity Detail Protein A affinity purified.

Species Reactivity Mouse

Specificity Recongizes a core functional region of mouse IL-4, which

prevents this cytokine from interacting with its cognate

receptor.

UniProt ID P07750

Last modified: May 29, 2024



info-

eu@enzolifesciences.com