

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 4 β -phorbol 12 β -myristate 12 α -acetate for 48h.
Isotype	IgG1 κ
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



Worry-free Guarantee

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +32 3 466 0420
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

This antibody is covered by our Worry-Free Guarantee