

IL-4 Recombinant monoclonal antibody (11B11) (Rabbit IgGκ)

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

ENZ-ABS548-0200	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	Rabbit
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.
Isotype	IgGκ
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.

Technical Info / Product Notes

This chimeric rabbit antibody was made using the variable domain sequences of the original Rat IgG1 format, for improved compatibility with existing reagents, assays and techniques.

UniProt ID

P07750

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-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