

ITAC monoclonal antibody (10)

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

Ordering Information

[Order Online »](#)

ENZ-ABS965-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Protected from light. Avoid freeze-thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CXCL11, H174, I-TAC, IP-9, C-X-C motif chemokine 11, Small-inducible cytokine B11, Interferon-inducible T-cell α chemoattractant, Interferon- γ -inducible protein 9, Beta-R1
Application	ELISA
Clone	10
Formulation	Liquid. 0.2 μ m filtered solution in PBS. Preservative free.
Host	Mouse
Immunogen	Recombinant human I-TAC / CXCL11 protein.
Isotype	IgG1
Purity Detail	Protein A purified.
Recommendation Dilutions/Conditions	ELISA (Capture) (1:250-1:2'000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES,
INC.
Phone: 800 942 0430
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(EU) 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

Life Sciences. September 17, 2024