

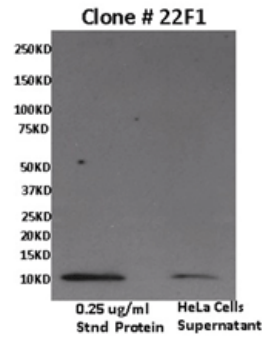
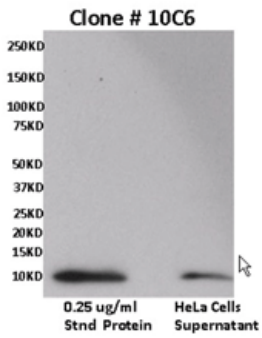
# ITAC monoclonal antibody (22F1) (biotin conjugate)

Interferon-inducible T-cell  $\alpha$  chemoattractant (ITAC), belonging to the CXC chemokine family, is expressed in peripheral blood leukocytes, pancreas and liver, thymus, spleen, lung, small intestine, placenta and prostate. Its molecular weight is 8–11 kDa. The cytokine is involved in regulating calcium release and inducing a chemotactic response in activated T-cells and is the dominant ligand for CXC receptor-3. The gene encoding this protein contains 4 exons and at least three polyadenylation signals which might reflect cell-specific regulation of expression. IFN- $\gamma$  is a potent inducer of transcription of this gene. It is upregulated in patients with AIDS, dementia and in patients with multiple sclerosis.

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

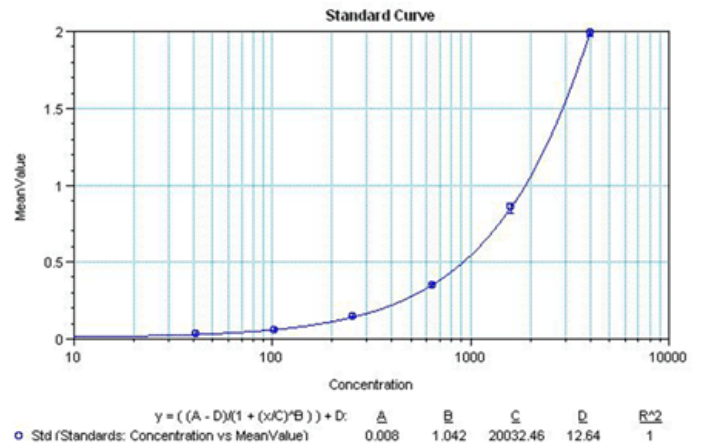
**Manuals, SDS & CofA**

[View Online »](#)



**Figure 1. Western Blot with ITAC, mAb (22F1) (biotin conjugate) (Prod no. ENZ-ABS275)**

Western blot analysis of human ITAC was performed by loading 0.25 µg of recombinant human ITAC and induced HeLa cell culture supernatant onto a 4-20% Tris-HCl polyacrylamide gel. Proteins were transferred to a PVDF membrane and blocked with 5% Milk/TBST for at least 1 hour. Membranes were probed with ITAC, mAb (22F1) (biotin conjugate) at a dilution of 1:1000 overnight at 4°C on a rocking platform. Membranes were washed in TBS-0.1% Tween 20 and probed with a goat anti-mouse-HRP secondary antibody. Membranes were washed and chemiluminescent detection performed.



**Figure 2. ELISA with ITAC, mAb (22F1) (biotin conjugate) (Prod no. ENZ-ABS275)**

Sandwich ELISA analysis of Human ITAC.

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	CXCL11, H174, I-TAC, IP-9, C-X-C motif chemokine 11, Small-inducible cytokine B11, Interferon-inducible T-cell $\alpha$ chemoattractant, Interferon- $\gamma$ -inducible protein 9, Beta-R1
<b>Application</b>	ELISA, WB
<b>Clone</b>	22F1
<b>Formulation</b>	Liquid. In PBS containing 4% BSA and a proprietary preservative.
<b>Host</b>	Mouse
<b>Immunogen</b>	Full length recombinant ITAC produced in E. coli.
<b>Isotype</b>	IgG1
<b>Purity Detail</b>	Protein G affinity purified.
<b>Recommendation Dilutions/Conditions</b>	ELISA (1 $\mu$ g/ml for coating and 0.125 – 0.25 $\mu$ g/ml for detection.)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	O14625
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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