

IFNAR1 polyclonal antibody

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

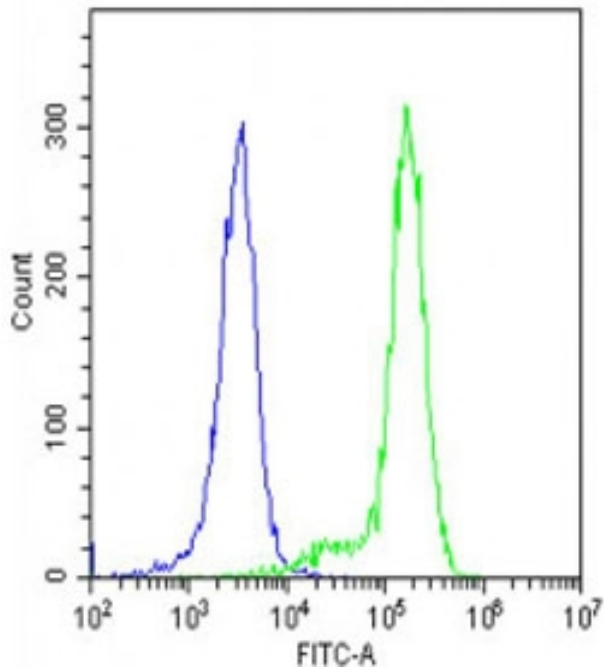
Ordering Information

[Order Online »](#)

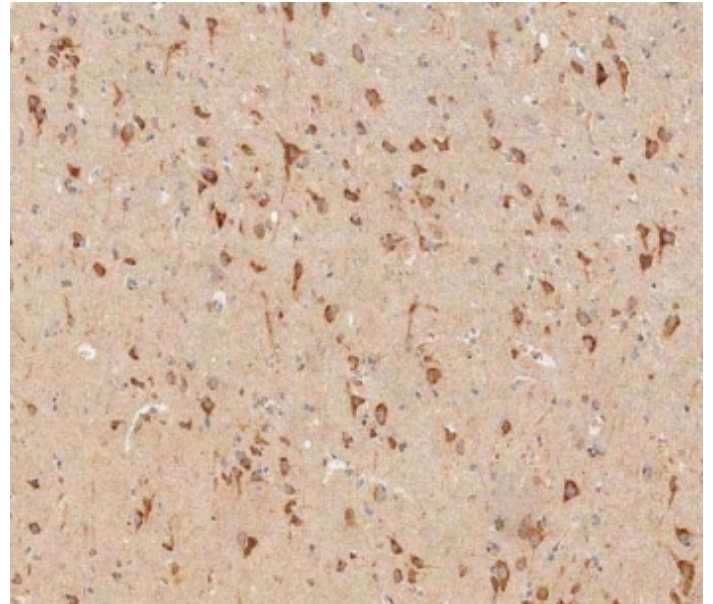
ENZ-ABS894-0400	0.4ml
-----------------	-------

Manuals, SDS & CofA

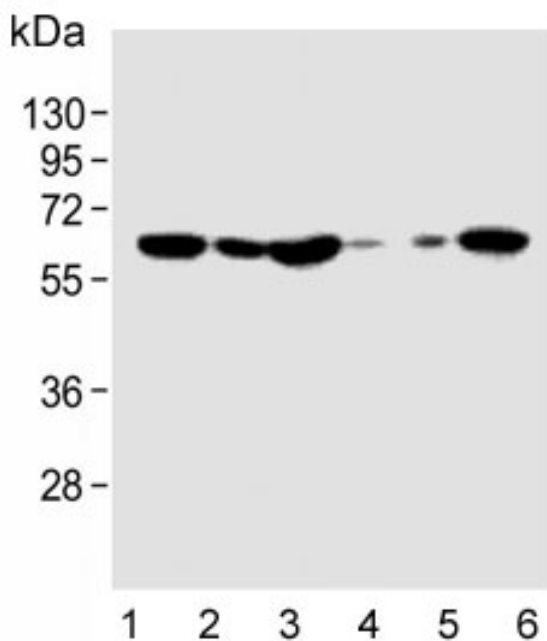
[View Online »](#)



Flow cytometry testing of fixed and permeabilized human K562 cells with Interferon alpha receptor 1 antibody; Blue=isotype control, Green= Interferon alpha receptor 1 antibody.



IHC testing of FFPE human brain tissue with Interferon alpha receptor 1 antibody. HIER: steam section in pH9 EDTA buffer for 20 min and allow to cool prior to staining.



Western blot testing of human 1) HeLa, 2) 293T, 3) Jurkat, 4) brain, 5) heart and 6) K562 cell lysate with Interferon alpha receptor 1 antibody. Expected molecular weight: 64-135 kDa depending on glycosylation level.

Handling & Storage

Handling	Aliquot the Interferon alpha receptor 1 antibody and store frozen at -20°C or colder. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Interferon- α/β receptor 1
Application	Flow Cytometry, IHC (PS), WB
Formulation	Liquid. In 1X PBS, pH 7.4, containing 0.09% sodium azide.
Host	Rabbit
Immunogen	Human IFNAR1 (aa 162-188).
Isotype	IgG
Purity Detail	Antigen affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (1:25, 1×10 ⁶ cells)Immunohistochemistry (FFPE) (1:50-1:100)Western blot (1:500-1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P17181
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



Last modified: July 12, 2024

ENZO LIFE SCIENCES,
INC.
Phone: +1 949 241 1200
usa@enzolifesciences.com

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

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