

# CD64 monoclonal antibody (AT37F7)

CD64 belongs to the immunoglobulin superfamily and it is a high-affinity Fc-gamma receptor. It plays a role in both innate and adaptive immune responses.

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

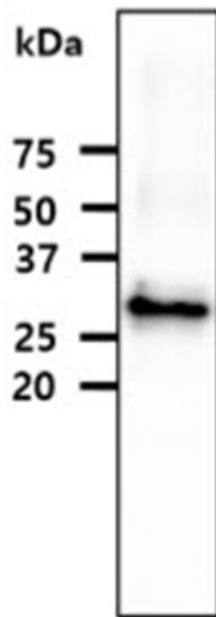
## Ordering Information

[Order Online »](#)

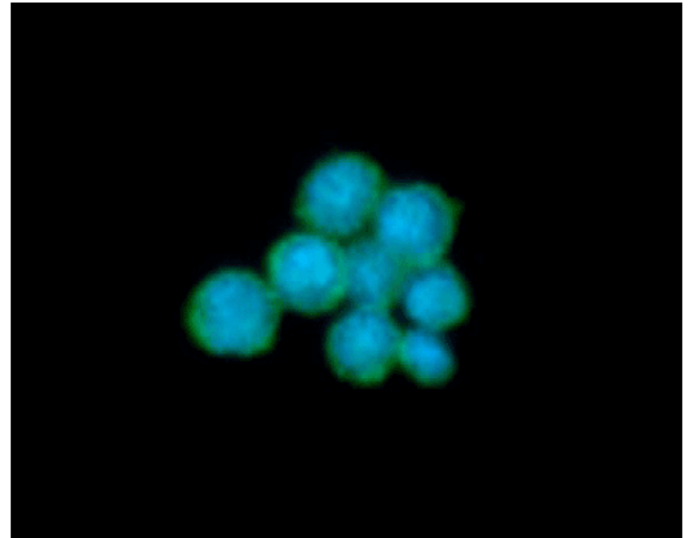
ENZ-ABS725-0100	100µl
-----------------	-------

## Manuals, SDS & CofA

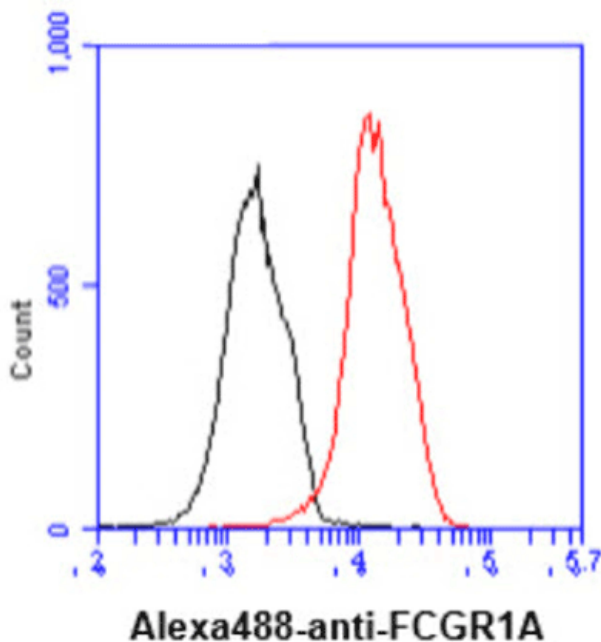
[View Online »](#)



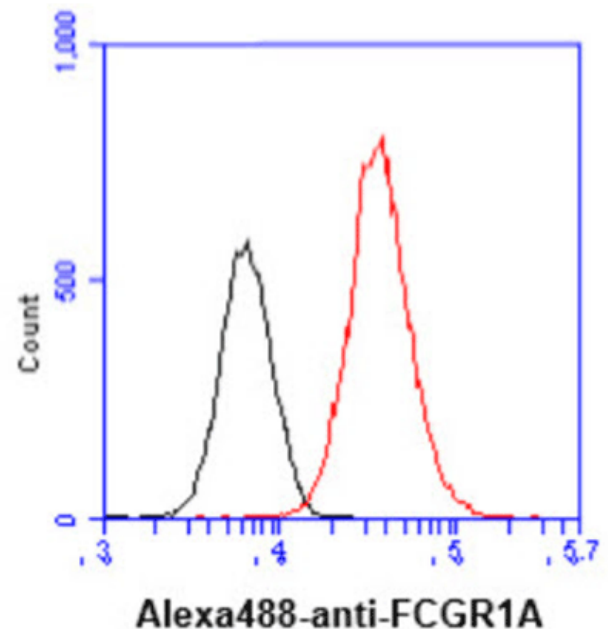
The recombinant protein (20µg) were resolved by SDS-PAGE, transferred to PVDF membrane and probed with anti-human CD64 antibody (1:1000). Proteins were visualized using a goat anti-mouse secondary antibody conjugated to HRP and an ECL detection system.



ICC/IF analysis of CD64 in Jurkat cells line, stained with DAPI (Blue) for nucleus staining and monoclonal anti-human FCGR1A antibody (1:100) with goat anti-mouse IgG-Alexa fluor 488 conjugate (Green).



Flow cytometry analysis of ENZ-ABS725 in Jurkat cell line, staining at 2-5µg for  $1 \times 10^6$  cells (red line). The secondary antibody used goat anti-mouse IgG Alexa fluor 488 conjugate. Isotype control antibody was mouse IgG (black line).



Flow cytometry analysis of ENZ-ABS725 in HeLa cell line, staining at 2-5µg for  $1 \times 10^6$  cells (red line). The secondary antibody used goat anti-mouse IgG Alexa fluor 488 conjugate. Isotype control antibody was mouse IgG (black line).

# Handling & Storage

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

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	High affinity immunoglobulin γ Fc receptor I, IgG Fc Receptor I, IGFR1, FcγRIa
Application	ELISA, Flow Cytometry, ICC, IF, WB
Clone	AT37F7
Formulation	Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.
Host	Mouse
Immunogen	Recombinant human FCGR1A (aa 16-292) purified from E. coli.
Isotype	IgG2bk
Purity Detail	Protein A affinity purified
Recommendation Dilutions/Conditions	Western Blot (1:2,000).Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from ascites.
Species Reactivity	Human
UniProt ID	P12314
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>



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)