

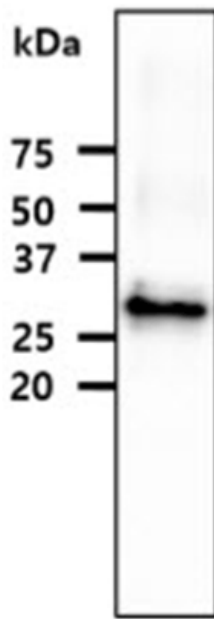
# CD64 monoclonal antibody (AT37F7)

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.

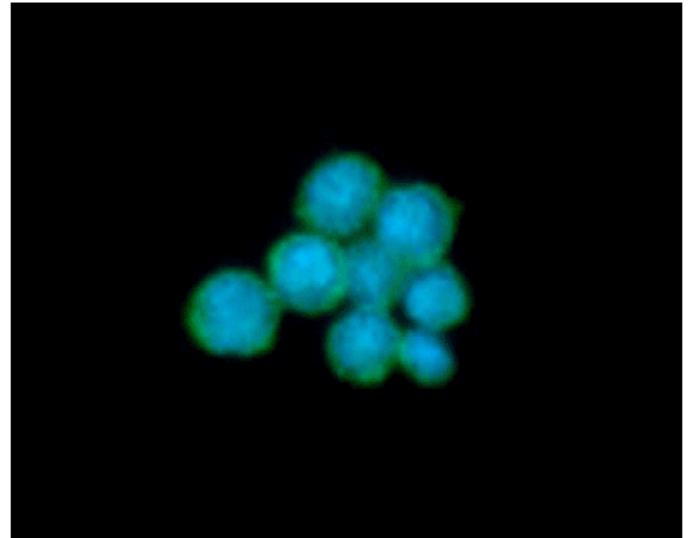
This antibody is covered by our [Worry-Free Guarantee](#).

**Manuals, SDS & CofA**

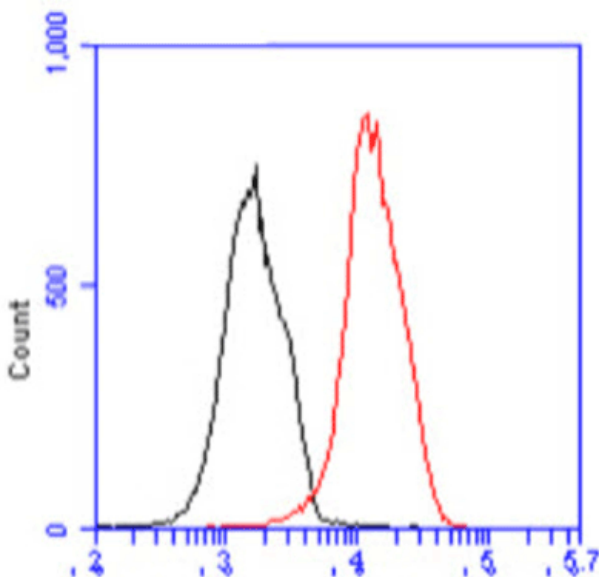
[View Online »](#)



The recombinant protein (20 $\mu$ g) 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.

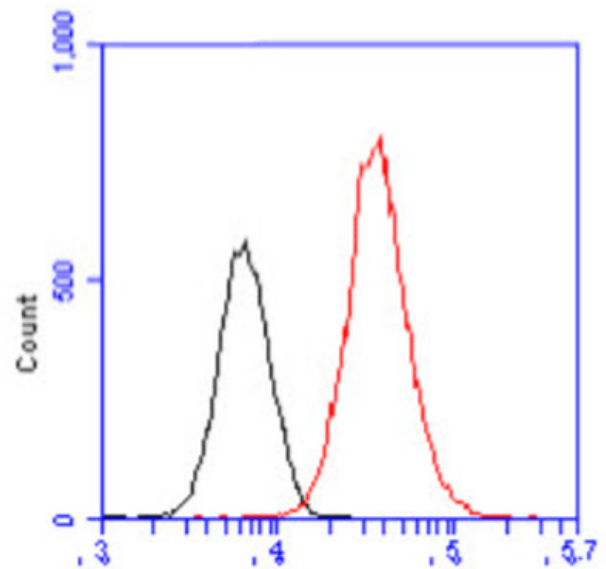


ICC/IF analysis of CD64 in Jurkat cells line, stained with DAPI (Blue) for nucleus staining and monoclonal anti-human FCGR1A antibody (1:100) with goat anti-mouse IgG-Alexa fluor 488 conjugate (Green).



**Alexa488-anti-FCGR1A**

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



**Alexa488-anti-FCGR1A**

Flow cytometry analysis of ENZ-ABS725 in HeLa cell line, staining at 2-5 $\mu$ g for 1 $\times$ 10<sup>6</sup> 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

<b>Alternative Name</b>	High affinity immunoglobulin $\gamma$ Fc receptor I, IgG Fc Receptor I, IGFR1, Fc $\gamma$ R1a
<b>Application</b>	ELISA, Flow Cytometry, ICC, IF, WB
<b>Clone</b>	AT37F7
<b>Formulation</b>	Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant human FCGR1A (aa 16-292) purified from E. coli.
<b>Isotype</b>	IgG2bk
<b>Purity Detail</b>	Protein A affinity purified
<b>Recommendation Dilutions/Conditions</b>	Western Blot (1:2,000). Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
<b>Source</b>	Purified from ascites.
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	P12314
<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)