

# RGMC (human) monoclonal antibody (1D7)

The RGMC (also called Hemojuvelin) is a GPI-anchored protein involved in iron metabolism through regulation of hepcidin expression. Mutations in the gene encoding RGMC cause severe iron overload (juvenile hemochromatosis /hemochromatosis type 2A). The protein binds to neogenin and acts as co-receptor of the bone morphogenetic protein. The mature RGMc protein is proteolytically processed and glycosylated.

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

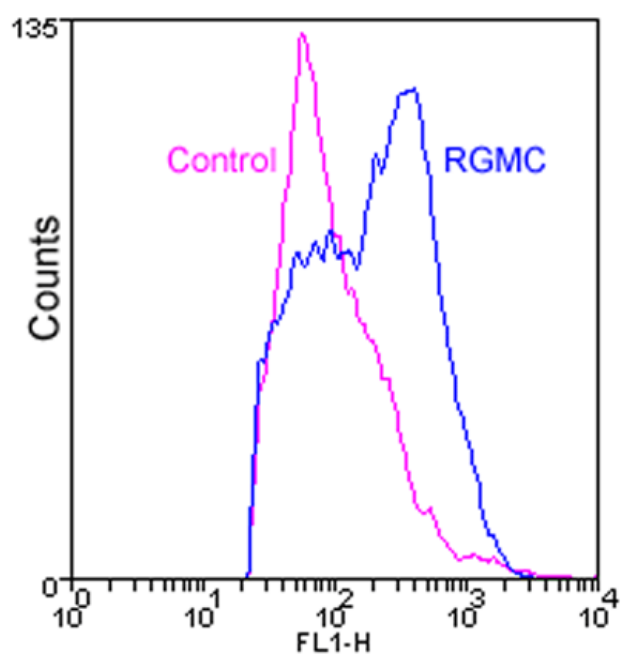
## Ordering Information

[Order Online »](#)

ENZ-ABS258-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of HEK293 cells stably overexpressing RGMC stained with RGMC (human), mAb (1D7) [Prod no. ENZ-ABS258] at 1:100 dilution (blue line) or normal mouse IgG (control, purple line).

# Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Repulsive guidance molecule C, Hemojuvelin, Hemochromatosis type 2 protein, HFE2, HJV
Application	ELISA, Flow Cytometry, IP, WB
Clone	1D7
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human RGMC (aa 36-398).
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Source	Purified from ascites.
Species Reactivity	Human
UniProt ID	Q6ZVN8
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)