

CD42a/CD42b/CD42c/CD42d complex (human) monoclonal antibody (Gi10)

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-805-023-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	GP1b/IX/V complex
Application	Flow Cytometry, ICC, IP
Clone	Gi10
Formulation	Liquid. 100µg of purified antibody in 200µl PBS containing 0.05% sodium azide.
Host	Mouse
Immunogen	Human platelets.
Isotype	IgG1
Species Reactivity	Human
Specificity	Recognizes the remnant moiety of the CD42a/CD42b/CD42c/CD42d (GP1b/IX/V) complex.
UniProt ID	P14770 (CD42a), P07359 (CD42b), P13224 (CD42c), P40197 (CD42c)
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .

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-be@enzolifesciences.com)

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