

CD42c (human) monoclonal antibody (Gi27)

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-805-026-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 β , GPIb β

Application IP, WB

Clone Gi27

Formulation Liquid. 100 μ g of purified antibody in 200 μ l PBS containing 0.05% sodium azide.

Host Mouse

Immunogen Human platelets.

Isotype IgG2b

Recommendation Dilutions/Conditions Western Blot (1:20)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

Specificity Recognizes the GPIb β subunit (CD42c) of the GPIb/IX/V (CD42a/CD42b/CD42c/CD42d) complex.

UniProt ID P13224

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

.