## CD49b/CD29 complex (human) monoclonal antibody (Gi14)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ALX-805-030-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 GPIa/lla complex

Application Antigen capture assay, Flow Cytometry, IHC (FS), IHC

(PS), IP

Clone Gi14

Formulation Liquid. Purified antibody in PBS containing 0.05% sodium

azide.

**Host** Mouse

**Immunogen** Human platelets.

lsotype lgG1

Species Reactivity Human

Specificity Recognizes the GPIa subunit (CD49b) of the GPIa/IIa

(CD49b/CD29) complex.

**UniProt ID** P17301 (CD49b), P05556 (CD29)

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

