G protein α_{i1} subunit monoclonal antibody (R4.5)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-SA281-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Gai1

Application WB

Clone R4.5

Crossreactivity Does not cross-react with transducin and subunits $G\alpha 0$,

Gai2, Gai3 or Gas.

Formulation Liquid. In PBS containing 0.2% BSA and 0.1% sodium

azide.

Host Mouse

Immunogen Partially purified rat brain Gαi1 protein.

Isotype IgG2b

Purity Detail Purified.

Species ReactivityBovine, Guinea pig, Human, Mouse, Rat

Specificity Recognizes the Ga_{i1} subunit of G protein.

UniProt ID P10824

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com

ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989

om infoeu@enzolifesciences.com

European Sales Office

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com

Last modified: May 29, 2024

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners