

# G protein $\alpha_{i1}$ 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 Ga0, Gai2, Gai3 or Gas.

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

**Host** Mouse

**Immunogen** Partially purified rat brain Gai1 protein.

**Isotype** IgG2b

**Purity Detail** Purified.

**Species Reactivity** Bovine, Guinea pig, Human, Mouse, Rat

**Specificity** Recognizes the  $G\alpha_{i1}$  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  
[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)