

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

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

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 Gα0, 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 Gα ₁₁ 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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Last modified: May 29, 2024