

G protein α_{i3} subunit polyclonal antibody

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Gai3
Application	IP, WB
Formulation	Liquid. In 100mM potassium phosphate, pH 7.5, containing 140mM sodium chloride. Contains no preservatives.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 345-354 of human Gai3 subunit.
Purity Detail	Protein A affinity purified IgG
Species Reactivity	Bovine, Human
Specificity	Recognizes the α_{i3} subunit of G protein.
UniProt ID	P08754
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



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-fr@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