

IgE (human) monoclonal antibody (I27) (Alexa Fluor[®] 647)

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

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage -20°C

Long Term Storage -80°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	Flow Cytometry
Clone	I27
Crossreactivity	Does not cross-react with human IgG, IgA or IgM.
Formulation	Liquid. In PBS containing 50% glycerol.
Host	Mouse
Immunogen	Purified human IgE (preincubated with mouse serum anti human IgG).
Isotype	IgG1
Species Reactivity	Human
Specificity	Recognizes human IgE (heat labile epitope in N-terminal of CH3 domain).
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-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