## IgE (human) monoclonal antibody (I27) (Alexa Fluor<sup>®</sup> 488)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

**ENZ-ABS213-0100** 

100µg

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 127

**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)

lsotype lgG1

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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

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

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com