Rat IgG2a Isotype Control (ATTO 488 conjugate)

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

Order Online »

ALX-804-836TD-C050

ATTO 488 - 50µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stable for at least

1 year after receipt when stored as recommended.

Handling Do not freeze. Protect from light.

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry

Application Notes Negative control in assays where Rat IgG (ATTO 488)-

labelled antibodies are used.

Clone KLH/G2a-11

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Rat

lgG2a κ

Species Reactivity Species independent

Technical Info / Product Notes ATTO-fluorescent antibodies show increased

photostability, outstanding brightness and intense signals. ATTO dyes are thermally stable, resistant to environmental changes and show no significant isomerization. ATTO 488 shows bright green fluorescence (λabs (max): 501nm, λem

This antibody is covered by our Worry-Free Guarantee

(max): 523nm, Emax: 90'000).

Worry-free Guarantee



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

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com

Last modified: May 29, 2024