

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
Isotype	IgG2a 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 (λ_{abs} (max): 501nm, λ_{em} (max): 523nm, ϵ_{max} : 90'000).

Worry-free Guarantee

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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

Last modified: May 29, 2024

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.