

Rabbit IgG Isotype control polyclonal antibody (DyLight488™ conjugate)

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	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 Rabbit IgG (DyLight™ 488)-labeled antibodies are used.
Emission Maximum	518 nm
Excitation Maximum	493 nm
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Host	Rabbit
Isotype	IgG
Source	Purified from rabbit serum.
Species Reactivity	Species independent
Technical Info / Product Notes	DyLight™ 488 has high fluorescence intensity over a broad pH range (pH 4-9) and is more photostable than many other dyes. The high water solubility of DyLight allows a high dye-to-protein ratio to be achieved without causing precipitation of the conjugates. Bright green fluorescence (λ_{abs} (max): 493nm, λ_{em} (max): 518nm, ϵ_{max} : 70,000).

Last modified: May 29, 2024



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-
be@enzolifesciences.com](mailto:info-be@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)