

Donkey anti-Rabbit IgG, Polyclonal Antibody (SEEBRIGHT® Red

594)

Donkey anti-Rabbit IgG polyclonal secondary antibody conjugated to SEEBRIGHT® Red 594 fluorophore, for detection of primary antibodies raised in rabbit in immunofluorescence (IF), immunocytochemistry (ICC), and immunohistochemistry (IHC) applications, including as a detection reagent option in Enzo fluorescence-based ISH/IHC workflows.

Fluorescently labeled secondary antibodies enable indirect immunodetection of unconjugated primary antibodies, providing signal amplification and multiplexing flexibility when paired with spectrally distinct fluorophores on other secondary antibodies.

Ordering Information

[Order Online »](#)

ENZ-ABS1075-0500	500µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

- Polyclonal donkey anti-Rabbit IgG, conjugated to SEEBRIGHT® Red 594
- Bright, photostable fluorescent signal for IF/ICC/IHC
- Minimal cross-reactivity for multiplex immunofluorescence panels

Handling & Storage

Use/Stability	Stable under recommended storage conditions up to the expiration date indicated on the vial label. Protect from light. Avoid repeated freeze-thaw cycles.
Handling	Protect from light. Centrifuge briefly before use. Avoid freeze-thaw cycles.
Short Term Storage	2–8°C
Long Term Storage	2–8°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	IF, ISH/IHC multiplex
Emission Maximum	617 nm
Excitation Maximum	590 nm
Host	Donkey
Recommendation Dilutions/Conditions	Optimal working dilution should be determined by the end user for each application.
Species Reactivity	Rabbit

Technical Info / Product Notes	Store at 2–8°C. Do not freeze. Centrifuge briefly before use. Optimal working dilution should be determined by titration for each specific application and tissue type. Compatible with AMPIVIEW® RNA-ISH workflows using SEEBRIGHT® fluorescence-based detection as an alternative to chromogenic (POLYVIEW® PLUS / HIGHDEF®) detection.
---	--



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)