

5/6-

Carboxyrhodamine 110-PEG3-Azide (25 mg)

With excitation maximum at 501 nm, the azide - carboxyrhodamine 110 conjugate is an excellent match to the intense 488 nm spectral line of the argon ion laser commonly used in many confocal laser-scanning microscopes. Carboxyrhodamine 110-PEG3-azide is an excellent green fluorescent rhodamine dye. The conjugate has better solubility than the parent dye carboxyrhodamine and its long PEG spacer should reduce any steric effect of the dye on the molecule to be labelled. Carboxyrhodamine 110 and its derivatives are, in general, significantly more photostable than any other known green fluorescent dyes including Alexa 488. In addition, the fluorescence of the azide-PEG3-carboxyrhodamine 110 conjugate is completely insensitive to pH between 4 and 9. Unlike Alexa 488, which is readily degraded under alkaline condition, carboxyrhodamine 110 and its derivatives are highly stable under both acidic and basic conditions. As a result, carboxyrhodamine 110 and its derivatives are superior alternatives to other green fluorescent dyes.

Ordering Information

[Order Online »](#)

JBS-CLK-AZ105-25

25mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance dark red solid

Couple Type Dye

Emission Maximum 525 nm

Excitation Maximum 501 nm

Formula $C_{29}H_{30}N_6O_7$

MW 574.59

Purity > 90 % (HPLC)

Solubility DMF, DMSO, MeOH

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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