

# 5/6-

# Carboxyrhodamine 110-PEG3-Azide (5 x 1 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-AZ105-5	5x1mg
-----------------	-------

## 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](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)