

5(6)-TRITC (ultra pure)

Amine reactive dye

5(6)-TRITC is an amino-reactive labeling reagent that is widely used in preparing bioconjugates of proteins and nucleic acids. The resultant conjugates have similar spectral properties to those prepared from 5(6)-TAMRA, SE with improved stability. TRITC is widely used and has a multitude of applications beyond protein, peptide and nucleotide labeling. For example, it can serve as an acceptor for FRET (Fluorescence Resonance Energy Transfer) based detection systems. FITC is recommended as a donor molecule in these FRET-based applications. Wavelength Maxima: Excitation 543nm, Emission 571nm

Citations: 3

[View Online »](#)

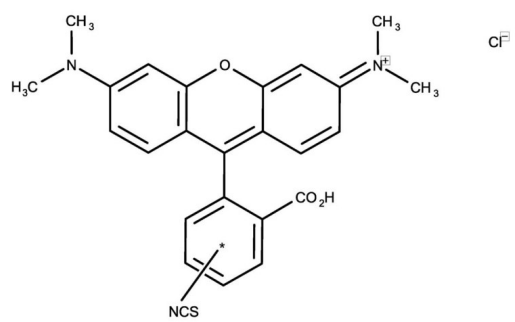
Ordering Information

[Order Online »](#)

ENZ-52452	10mg
-----------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended.
Handling	Protect from light. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Tetramethylrhodamine-5(6)-isothiocyanate
CAS	95197-95-8
Formula	$C_{25}H_{21}N_3O_3S$
MW	480.0
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.

Technical Info / Product Notes	This product is a member of the CELLESTIAL [®] product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL [®] reagents and kits are optimal for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required.
--------------------------------	---



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)