

5-FITC (ultra pure)

Amine reactive dye

Fluorescein isothiocyanate (FITC) is an amine-reactive fluorescein derivative that yields conjugates with superior stability. Dye synthesis generally results in two isomers that differ slightly in protein binding and may elute differently under chromatographic conditions, complicating analysis in certain situations. The 5-isomer of FITC is the most widely used isoform and has a multitude of applications beyond protein peptide and nucleotide labeling. For example, it is commonly employed as a donor for FRET (Fluorescence Resonance Energy Transfer) based assay systems. An acceptor molecule, such as TRITC, is recommended in such FRET applications. Wavelength Maxima: Excitation 494nm, Emission 520nm

Citations: 2

[View Online »](#)

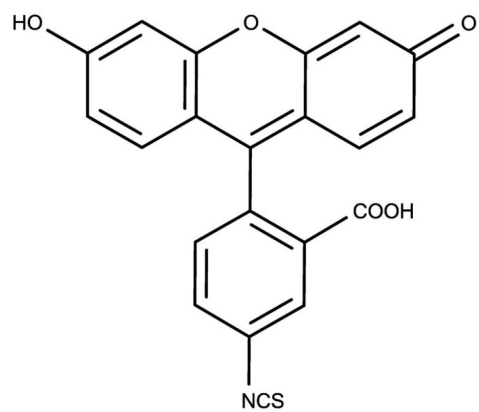
Ordering Information

[Order Online »](#)

ENZ-52451	1g
-----------	----

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	5-Fluorescein Isothiocyanate
CAS	3326-32-7
Formula	$C_{21}H_{11}NO_5S$
MW	389.4
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)