

Fluorescein-5-maleimide (ultra pure)

Thiol dye

Fluorescein-5-maleimide enables protein labeling via thiol modifications in most proteins at cysteine residues that either are intrinsically present or resulted from reduction of cystine residues. Unlike iodoacetamides, maleimides do not react with histidine and methionine under physiological conditions. The probe is suitable for assessing conformation- dependent accessibility of cysteine residues in proteins. Wavelength Maxima: Excitation 493nm, Emission 515nm

Citations: 2

[View Online »](#)

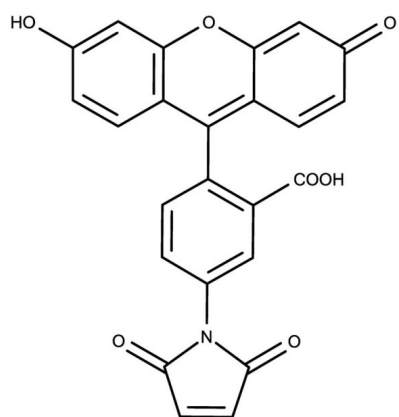
Ordering Information

[Order Online »](#)

ENZ-52502	25mg
-----------	------

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

Appearance	Orange solid.
CAS	75350-46-8
Formula	$C_{24}H_{13}NO_7$
MW	427.4
Purity	≥90% (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)