

# Nile Red (ultra pure)

## Lipid dye

Nile red is a lipophilic probe that becomes intensely fluorescent in a hydrophobic environment, while it has minimal fluorescence in aqueous media. It is an excellent vital stain for the detection of intracellular lipid droplets by fluorescence microscopy, in cells such as adipocytes. Flow cytometry methods using a single staining with Nile red and double staining with Nile red and anti-CD3 monoclonal antibody have been employed to monitor peripheral leukocyte and lymphocyte phospholipidosis. Nile Red has also been used for rapid staining of proteins in SDS-polyacrylamide gel electrophoresis. Wavelength Maxima: Excitation 552nm, Emission 636nm

Citations: 3

[View Online »](#)

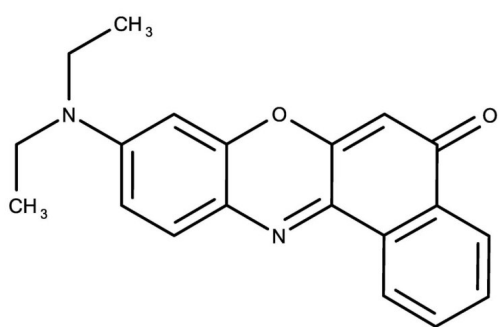
## Ordering Information

[Order Online »](#)

ENZ-52551	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

CAS	7385-67-3
MW	318.4
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.

## Technical Info / Product Notes

This product is a member of the CELLESTIAL<sup>®</sup> product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL<sup>®</sup> 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.

Last modified: May 29, 2024



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)