

Dihydrorhodamine 123 (ultra pure)

Mitochondria dye

Dihydrorhodamine 123 is a widely utilized probe for detection of intracellular reactive oxygen species (ROS) such as peroxide, hypochlorous acid and peroxynitrite. It is readily oxidized into rhodamine 123, which exhibits a spectral profile similar to that of FITC. In combination with other fluorescent probes (such as surface receptor-targeted fluorescent antibodies, the cell viability probe propidium iodide, or fluorescent calcium indicators) dihydrorhodamine 123 can be used for multiparametric cell measurements. Wavelength Maxima: Excitation 507nm, Emission 529nm

Citations: 3

[View Online »](#)

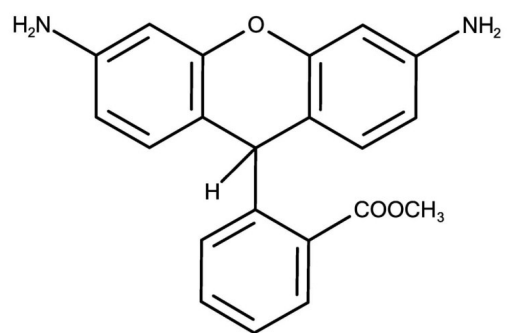
Ordering Information

[Order Online »](#)

ENZ-52302	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	2-(3,6-Diamino-9H-xanthene-9-yl)-benzoic acid methyl ester
CAS	109244-58-8
MW	346.4
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.
Technical Info / Product Notes	<p>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.</p>

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com