

Rhodamine 123 (ultra pure)

Mitochondria dye

Rhodamine 123 is a cell-permeant, cationic, green-fluorescent dye that is readily sequestered by active mitochondria without cytotoxic effects. Rhodamine 123 can be used in multi-parametric analysis, without fluorescence interference, in combination with common protein labeling dyes such as cyanine-5 and AMCA. With respect to apoptosis, the presence of mitochondrial membrane potential can be probed with rhodamine 123 while the structure and integrity of mitochondria can be assessed using 10-N-nonyl-acridine orange. Wavelength Maxima: Excitation 507nm, Emission 529nm

Citations: 8

[View Online »](#)

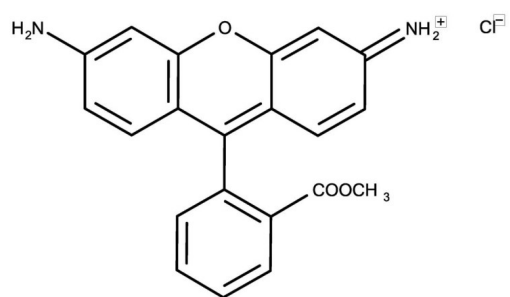
Ordering Information

[Order Online »](#)

ENZ-52307	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

Alternative Name	3,6-Diamino-9-(2-(methoxycarbonyl)phenyl)xanthylium chloride
CAS	62669-70-9
Formula	$C_{21}H_{17}ClN_2O_3$
MW	380.8
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-fr@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