

TMRE (ultra pure)

Mitochondria dye

Positively charged rhodamine dyes (such as rhodamine esters and rosamines) are selectively localized in mitochondria, thus they are widely used for labeling mitochondria of live cells. Like JC-1, TMRE is widely used for measuring mitochondrial membrane potential, in addition to selectively staining mitochondria. Real-time imaging of mitochondrial membrane potential in individual cardiomyocytes within perfused rat hearts has been demonstrated with this dye, using 2-photon laser-scanning microscopy. Wavelength Maxima: Excitation 549nm, Emission 574nm

Citations: 15

[View Online »](#)

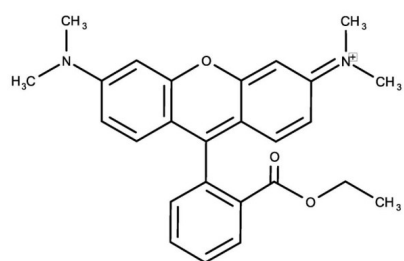
Ordering Information

[Order Online »](#)

ENZ-52309	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	Tetramethylrhodamine ethyl ester perchlorate, 3,6-bis(Dimethylamino)-9-[2-(ethoxycarbonyl)phenyl]-xanthylium perchlorate
CAS	115532-52-0
Formula	$C_{26}H_{27}ClN_2O_7$
MW	515.0
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

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