

# DiIC12(3) perchlorate (ultra pure)

## Membrane dye

DiI, DiO, DiD and DiR dyes are a family of lipophilic fluorescent stains for labeling membranes and other hydrophobic structures. The fluorescence of these environmentsensitive dyes is greatly enhanced when incorporated into membranes or bound to lipophilic biomolecules such as proteins although they are weakly fluorescent in water. Once applied to cells, these dyes diffuse laterally within the cellular plasma membranes, resulting in even staining of the entire cell at their optimal concentrations. DiIC12(3) is a lipophilic neuronal tracer that is commonly used for labeling neuronal projections as well as lipid bilayers in other cell types, particularly endothelial cells. DiIC12(3) exhibits low toxicity and minimally effects cell viability and is tolerable in culture for several days, making it suitable across a variety of applications. Wavelength Maxima: Excitation 549nm, Emission 565nm

Citations: 7

[View Online »](#)

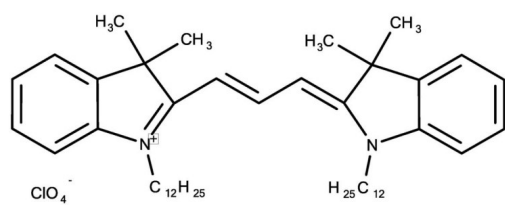
## Ordering Information

[Order Online »](#)

ENZ-52206	100mg
-----------	-------

## 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	1,1'-Didodecyl-3,3',3'-tetramethylindocarbocyanine perchlorate
Appearance	Dark red powder.
CAS	75664-01-6
Formula	$C_{47}H_{73}N_2O_4Cl$
MW	765.5
Purity	≥95% (TLC)
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.



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)