

# Propidium Iodide (ultra pure)

## DNA dye

Propidium iodide (PI) belongs to the same chemical class as ethidium bromide. As with ethidium bromide, its fluorescence is enhanced 20-30-fold upon binding to nucleic acids. The fluorescence excitation maximum is red-shifted 30-40 nm and the fluorescence emission maximum blueshifted ~15 nm. PI also binds to RNA as does DAPI and acridine orange. PI is membrane impermeant and is commonly used for identifying dead cells in a mixed population of cells and as a counter-stain in multicolor fluorescent techniques. It can also be used to differentiate necrotic, apoptotic and viable cells. The dye is suitable for fluorescence microscopy, flow cytometry and fluorometry applications. Wavelength Maxima: Excitation 535nm, Emission 617nm

Citations: 7

[View Online »](#)

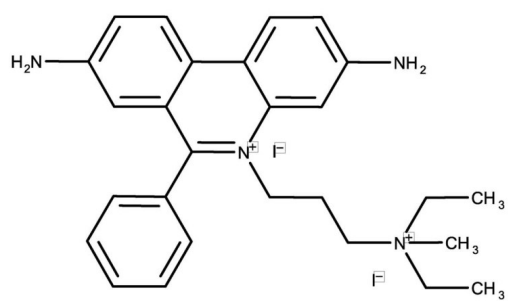
## Ordering Information

[Order Online »](#)

ENZ-52403	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

Appearance	Dark red solid.
CAS	25535-16-4
Formula	$C_{27}H_{34}I_2N_4$
MW	668.4
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.
--------------------------------	---



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)