

# DIBAC4(3) (ultra pure)

## Membrane Potential Detector

DiBAC4(3) is a sensitive slow-response probe for measuring cellular membrane potential. In general, slow-response probes exhibit potential-dependent changes in their transmembrane distribution that are accompanied by a fluorescence change. The magnitude of response is much larger than fast-response probes. Slow-response probes, which include cationic carbocyanines, rhodamines and anionic oxonols, are suitable for detecting changes in average membrane potentials of nonexcitable cells caused by respiratory activity, ion-channel permeability, drug binding and other factors. DiBAC4(3) has been employed in flow cytometry to monitor antibacterial activity of defensins. Wavelength Maxima: Excitation 493nm, Emission 516nm

Citations: 3

[View Online »](#)

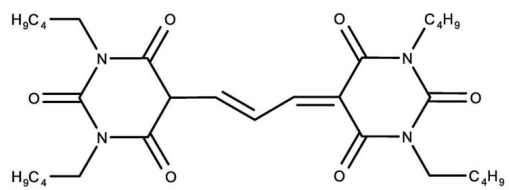
## Ordering Information

[Order Online »](#)

ENZ-52205	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	bis-(1,3-Dibarbituric acid)-trimethine oxonol
CAS	70363-83-6
Formula	$C_{27}H_{40}N_4O_6$
MW	516.6
Purity	≥98% (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)