

MQAE (ultra pure)

Chloride dye

MQAE is a fluorescent chloride ion indicator that is more sensitive and selective than ^{36}Cl and microelectrode-based methods for chloride measurement in cells. MQAE is a fluorescent indicator that is quenched via collision with chloride. The probe is suitable for detection of chloride concentrations in liposomes, reconstituted membranes or in extracellular medium. As a UV-excitabile indicator, MEQ can be detected using an argon-ion laser for confocal microscopy and flow cytometry applications. MQAE has greater sensitivity than SPQ and higher fluorescence quantum yield. The ester group of MQAE may slowly hydrolyze inside cells, resulting in a change in its fluorescence response. Wavelength Maxima: Excitation 350nm, Emission 460nm

Citations: 2

[View Online »](#)

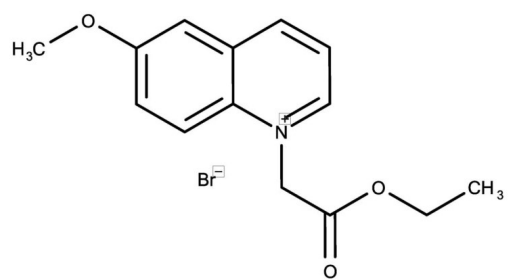
Ordering Information

[Order Online »](#)

ENZ-52156	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	N-(Ethoxycarbonylmethyl)-6-methoxyquinolinium bromide
CAS	162558-52-3
Formula	$C_{14}H_{16}BrNO_3$
MW	326.2
Purity	≥98% (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](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)