

TSQ (ultra pure)

Zinc dye

TSQ is an efficient membrane-permeable fluorescent probe for zinc (II). It is selective for Zn (II) in the presence of physiological concentrations of Ca (II) and Mg (II) ions. Fluorescence of TSQ is enhanced upon binding Zn (II), but the probe does not demonstrate a shift or change in emission wavelength. TSQ forms a largely insoluble 2:1 (Ligand-Metal) complex with zinc and emits blue light upon excitation at 365 nanometers.

Citations: 14

[View Online »](#)

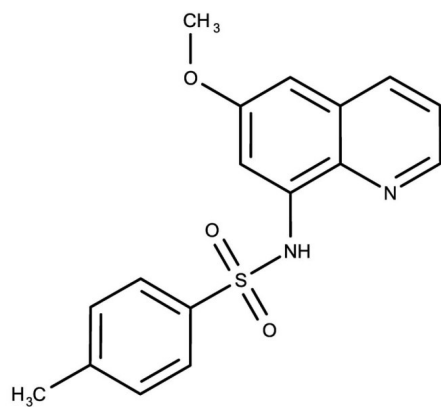
Ordering Information

[Order Online »](#)

ENZ-52153	25mg
-----------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended. In solvent: – 80°C, 6 months; -20°C, 1 month.
Handling	Protect from light. Keep cool and dry.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name N-(6-Methoxy-8-quinolyl)-p-toluenesulfonamide

Appearance Light brown to brown solid.

CAS 109628-27-5

Emission Maximum 385 nm

Excitation Maximum 344 nm

Formula $C_{17}H_{16}N_2O_3S$

MW 328.4

Purity ≥98% (LCMS)

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-
fr@enzolifesciences.com](mailto:info-fr@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)