

# 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 »](#)

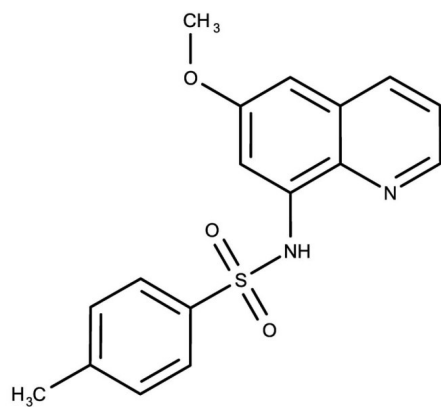
## Ordering Information

[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<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-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)

Last modified: November 5, 2024

For local distributors and detailed product information visit us online: [enzo.com](http://enzo.com) All product names, logos, and brands are the property of their respective owners.