SPQ

The fluorescence of SPQ is specifically quenched by chloride via collision. Therefore, chloride concentration is measured by monitoring the degree of fluorescence decrease.



Product attributes

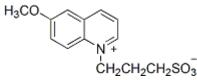
CAS number	83907-40-8
Cell permeability	Membrane impermeant
Colors	Blue
Excitation/Emission	344/443 nm

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

The fluorescence of SPQ is specifically quenched by chloride via collision. Therefore, chloride concentration is measured by monitoring the degree of fluorescence decrease. The dye can be loaded into cells by hypotonic shock.

- White solid soluble in H2O and DMSO
- Store at 4°C. Protect from light, especially in solution.
- \bullet $\lambda_{Ex}/\lambda_{Em}=344/443$ nm
- ϵ (344 nm, H₂O) = 3,700
- C₁₅H₁₅NO₄S
- MW: 281
- [83907-40-8]



As the indicator does not covalently bind to cellular components, it may be actively effluxed from the cell by organic anion transporters. The rate of efflux increases with temperature, and may vary between cell types, resulting in variable retention times of a few minutes to hours. Experiments using indicators in cells usually are performed within one or two hours of loading, but it may be possible to re-load cells with indicator if needed. The organic anion transporter inhibitor Probenecid (#50027) can be used to slow the rate of indicator efflux from cells.

References

- 1. Neurosci Lett, 104, 326 (1989), DOI: 10.1016/0304-3940(89)90597-1
- 2. Biophys J, 56, 1071 (1989), DOI: 10.1016/S0006-3495(89)82755-9
 3. J Biol Chem, 266, 20590 (1991), DOI: 10.1016/S0021-9258(18)54749-9
- 4. Methods Cell Biol, 99, 113, (2021), DOI: 10.1016/B978-0-12-374841-6.00005-0

This datasheet was generated on November 9, 2025 at 08:40:42 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/spq6-methoxy-n-3-sulfopropylquinolinium/