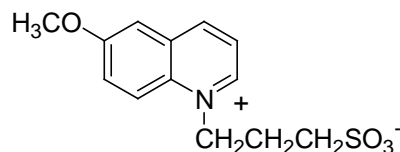


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **SPQ** (6-methoxy-N-(3-sulfopropyl)quinolinium)**CATALOG #:** **52010****MOLECULAR INFORMATION:** C₁₃H₁₅NO₄S
Mwt: 281
[83907-40-8]**PROPERTIES:**

Color & Form	White solid
Purity	≥95% by HPLC
Solubility	Soluble in water and DMSO
λ_{ex}/λ_{em}	344nm/443nm
ε	3,000

STORAGE AND HANDLING: Store at 4 °C. Protect material from light, especially in solution**APPLICATION:** 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.

Ref: 1) Engblom, A.C., et al. Neurosci. Lett. 104, 326(1989); 2) Chao, A.C., et al. Biophys. J. 56, 1071(1989); 3) Mulberg, A.E., et al. J. Biol. Chem. 266, 20590(1991)

TOXICITY: Unknown.

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--