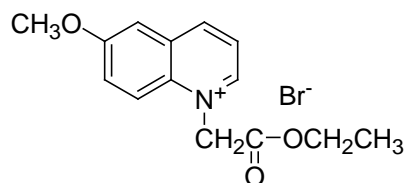


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: MQAE (N-(ethoxycarbonylmethyl)-6-methoxyquinolinium bromide)**CATALOG #:** 52011**MOLECULAR INFORMATION:** C₁₄H₁₆BrNO₃
Mwt: 326.19**PROPERTIES:**

Color & Form	Light brown solid
Purity	≥99% by HPLC
Solubility	Soluble in water (≥1 mg/mL)
$\lambda_{ex}/\lambda_{em}$	350nm/460nm
ϵ	2,800

STORAGE AND HANDLING: Store at 4 °C. Protect material from light, especially in solution**APPLICATION:** MQAE is an improved chloride indicator that has greater sensitivity to chloride ($K_{sv}=200 \text{ M}^{-1}$) than SPQ ($K_{sv}=118 \text{ M}^{-1}$) and higher fluorescence quantum yield. The ester group of MQAE may slowly hydrolyze inside cells, resulting in a change in its fluorescence response.Ref: 1) *Am J Physiol* **259**, C375 (1990); 2) *Anal Biochem* **178**, 355 (1989); 3) *Anal Biochem* **241**, 51(1996)**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>
--