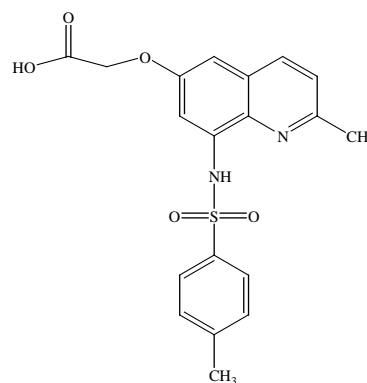


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Zinquin free acid**CATALOG #** 52022**MOLECULAR INFORMATION:** C₁₉H₁₈N₂O₅S
Mwt: 386.42**PROPERTIES:**

Color & Form	Off-white solid
Purity	≥ 98% by HPLC
Solubility	Soluble in DMSO
Absorption/Emission	364 nm/485 nm

STORAGE AND HANDLING: Stored at 4 °C upon receipt. Protect from light, especially when in solution.**APPLICATION:** Zinquin is an UV-excitable fluorescent zinc indicator. Zinc is believed to be involved in the suppression of apoptosis and play important roles in many neural activities.Ref.: 1) *Biochem. J.* **303**, 781(1994); 2) *Biochem. J.* **296**, 403(1993).**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>
--