

Instruction Manual

| | |
|---------------------------------------|---|
| Catalog Number | PK-CA707-52022 |
| Description | Zinquin is an UV-excitable, blue fluorescent zinc indicator. Zinc is believed to be involved in the suppression of apoptosis and play important roles in many neural activities. |
| Quantity | 5 mg |
| Excitation / Emission Maxima | $\lambda_{exc}/\lambda_{em}$ (in the presence of Zn^{2+}) = 364/385 nm |
| Molecular Structure | |
| Molecular Weight / Molecular Formula | 386.42 Da; $C_{19}H_{18}N_2O_5S$ |
| Purity | >98% (as determined by HPLC) |
| Appearance / Formulation / Solubility | Off-white solid; soluble in DMSO. |
| Storage & Stability | Store at 4°C. Protect from light, especially when in solution. |
| Applications | Fluorescent zinc indicator |
| References | 1) Biochem. J. 303, 781(1994) 2) Biochem. J. 296, 403(1993) |
| Caution | 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. |

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.