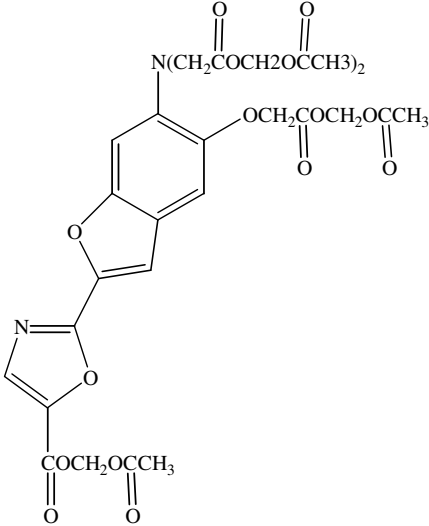


### Instruction Manual

|                                       |   |
|---------------------------------------|---|
| Catalog Number                        | PK-CA707-50037  |
| Description                           | Furaptra, AM ester is a membrane-permeant, UV light-excitable intracellular magnesium indicator (also known as Mag-Fura-2) and useful for noninvasive intracellular loading. The indicator is loaded into cells via simple incubation of cells in the buffer containing the dye. See information for Mag-Fura-2/tetrapotassium salt (PK-CA707-50035).           |
| Quantity                              | 10 x 100 µg   |
| Excitation / Emission Maxima          | $\lambda_{ex}/\lambda_{em} = 366/475$ nm (in EtOAc);<br>Extinction Coefficient: $31,000 \text{ M}^{-1}\text{cm}^{-1}$   |
| Molecular Structure                   |   |
| Molecular Weight / Molecular Formula  | 723 Da; $\text{C}_{30}\text{H}_{30}\text{N}_2\text{O}_{19}$   |
| Purity                                | >95% (as determined by HPLC)  |
| Appearance / Formulation / Solubility | Light yellow solid; soluble in DMSO.  |
| Storage & Stability                   | Store desiccated at $-20^\circ\text{C}$ . Protect from light.   |
| Applications                          | Fluorescent calcium and magnesium indicator   |
| References                            | <ol style="list-style-type: none"> <li>1) J Chromatogr A 1995; (696):93-93</li> <li>2) J Photochem Photobiol B 1996; (33):125-125</li> <li>3) Front Biosci 1997; (2):13</li> <li>4) Cell Biology: A Laboratory Handbook, 2nd Edn. (JE Celis, Ed.) 1998:363-374</li> <li>5) Cancer Cell 2002; (1):169-169</li> </ol>   |
| 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.