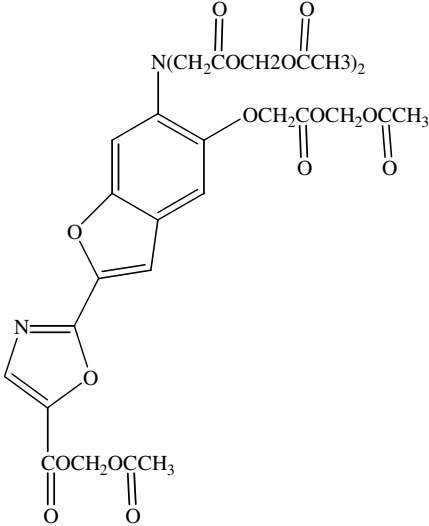


Instruction Manual

Catalog Number	PK-CA707-50039
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	20 x 50 µg
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em} = 366/475 \text{ nm}$ (in EtOAc); 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°C . Protect from light.
Applications	Fluorescent calcium and magnesium indicator
References	<ol style="list-style-type: none"> 1) J Chromatogr A 1995; (696):93-93 2) J Photochem Photobiol B 1996; (33):125-125 3) Front Biosci 1997; (2):13 4) Cell Biology: A Laboratory Handbook, 2nd Edn. (JE Celis, Ed.) 1998:363-374 5) Cancer Cell 2002; (1):169-169
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.