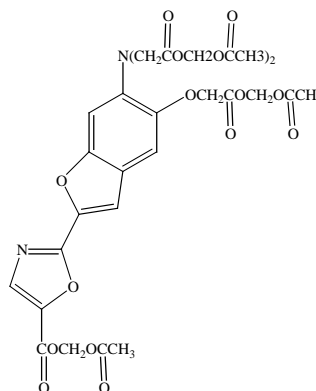


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **Furaptra, AM ester** (also known as **Mag-Fura-2, AM ester**)**CATALOG #:** **50037/50038/50039****MOLECULAR INFORMATION:** C₃₀H₃₀N₂O₁₉
MW: 723**PROPERTIES:**

Color & Form	Light yellow solid
Purity	≥ 95% by TLC
Solubility	Soluble in DMSO
Absorption/Emission	366nm/475nm (EtOAc)
Extinction Coefficient	31,000

STORAGE AND HANDLING: Store desiccated at -20 °C. Protect from light.**APPLICATION:** Membrane-permeant form of Furaptra. 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 (**50035**).**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>
--