

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

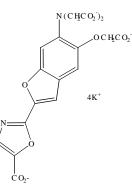
PRODUCT NAME:

Furaptra, tetrapotassium salt (also knowns as Mag-Fura-2, tetrapotassium salt)

CATALOG #:

MOLECULAR INFORMATION: $\begin{array}{c} C_{18}H_{10}K_4N_2O_{11} \\ MW: 587 \end{array}$

50035



PROPERTIES: Color & Form Light yellow solid Purity >95% by HPLC **Solubility** Soluble in water (pH > 6)**Absorption/Emission** 369nm/511nm (H₂O) Extinction 22,000 Coefficient STORAGE AND Store desiccated at -20 °C. Protect from light. HANDLING: Furaptra is a UV-excitable fluorescent indicator for magnesium with a K_d of 1.9 mM.^{1,2} **APPLICATION:** Similar to Fura-2, the excitation wavelength of Mag-Fura-2 undergoes a blue shift from 369 nm to 330 nm. Mag-Fura-2 also responds to Ca^{2+} but with a significantly lower K_d than Fura-2 to Ca^{2+} . An important application of Mag-Fura-2 is its use in detecting high, transient Ca^{2+} concentration during Ca^{2+} spikes.^{3,4} Mag-Fura-2/tetrapotassium salt can be loaded into cells by microinjection or scrape loading. Ref: 1)Pro. Natl. Acad. Sci. USA 86, 2981(1989); 2) Am. J. Physiol. 256, C540(1989); 3) Neuron 10, 21(1993); 4) Biophys. J. 68, 2156(1995). **TOXICITY:** Unknown. Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on FIRST AID: 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: *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.*