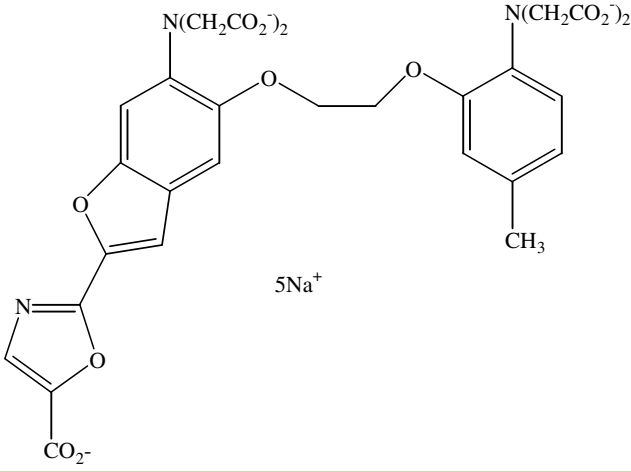


Fura-2 (pentasodium salt)

1-[6-Amino-2-(5-carboxy-2-oxazolyl)-5-benzofuranyloxy]-2-(2'-amino-5'-methylphenoxy)-ethane-N,N,N',N'-tetraacetic acid; pentasodium salt

Instruction Manual

Catalog Number	PK-CA707-50032
Description	Fura-2 is a widely used UV-excitable fluorescent calcium indicator developed by professor Roger Tsien. It has been used in many cellular systems and applications particularly in microscopic imaging. Upon calcium binding, the fluorescent excitation maximum of the indicator undergoes a blue shift from 363 nm (Ca ²⁺ -free) to 335 nm (Ca ²⁺ -saturated), while the fluorescence emission maximum is relatively unchanged at ~510 nm. The indicator is typically excited at 340 nm and 380 nm respectively and the ratio of the fluorescent intensities corresponding to the two excitations is used in calculating the intracellular concentrations. Measurement of calcium concentration using this ratioing method avoids interference due to uneven dye distribution and photobleaching. Fura-2 ammonium, potassium, or sodium salt is membrane-impermeant but can be loaded into cells via microinjection or scrape loading.
Quantity	1 mg
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em} = 363/512$ nm (no Ca ²⁺); $\lambda_{ex} \backslash \lambda_{em} = 335/505$ nm (high Ca ²⁺); Extinction Coefficient: 27,000 M ⁻¹ cm ⁻¹ (363 nm, no Ca ²⁺); 35,000 M ⁻¹ cm ⁻¹ (high Ca ²⁺)
Molecular Structure	 <p>5Na⁺</p>
Molecular Weight / Molecular Formula	752 Da; C ₂₉ H ₂₂ N ₃ Na ₅ O ₁₄
Purity	>95% (as determined by HPLC)
Appearance / Formulation / Solubility	Light yellow solid; soluble in water or DMSO.
Storage & Stability	Stored at ≤4°C upon receipt. Protect from light, especially when in solution.
Intended Use	For in vitro research use only. Not for diagnostic or therapeutic procedures.
Applications	Fluorescent calcium indicator The K _d for Fura-2 was reported to be 224 nM in cell-free media. However, the K _d is usually affected by a number of factors in cells including pH, proteins concentrations, ionic strength, temperature and viscosity. Thus, calibration of the K _d is necessary for accurate measurement of intracellular calcium concentrations. For details on calibration, we recommend that you consult the references listed at the end of this document (See refs 2-8). PromoKine offers A-23187 (PK-CA707-59001), an ionophore that is commonly used for intracellular calibration of calcium indicators. PromoKine also offers EDC (PK-CA707-59002, also known as EDAC), which can be used to fix calcium indicators in cells, if post histochemical studies are desired following physiological experiments.

References	<p>1) J. Biol. Chem. 260, 3440(1985) 2) Bright, G.R., et al, in Fluorescence Microscopy of Living Cells in Culture, Part B, (Methods in Cell Biology, Vol. 30), Academic Press (1989) p. 157 3) Am. J. Physiol. 261, C1107(1991) 4) Biophys. J. 54, 1089(1988) 5) Biochem. Biophys. Res. Comm. 177, 184(1991) 6) Cell Calcium 11, 85(1990) 7) Cell Calcium 12, 279(1991) 8) Neuropharmacol. 34, 1423(1995)</p>
Caution	<p>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.</p>

PromoCell GmbH

Sickingenstr. 63/65
69126 Heidelberg
Germany

Email: info@promokine.info
www.promokine.info

North America

Phone: 1 – 866 – 251 – 2860 (toll free)
Fax: 1 – 866 – 827 – 9219 (toll free)

Deutschland

Telefon: 0800 – 776 66 23 (gebührenfrei)
Fax: 0800 – 100 83 06 (gebührenfrei)

France

Téléphone: 0800 90 93 32 (ligne verte)
Téléfax: 0800 90 27 36 (ligne verte)

United Kingdom

Phone: 0800 – 96 03 33 (toll free)
Fax: 0800 – 169 85 54 (toll free)

Other Countries

Phone: +49 6221 – 649 34 0
Fax: +49 6221 – 649 34 40