

# PRODUCT INFORMATION



## Fura-2 (sodium salt)

Item No. 20414

**Formal Name:** 2,2'-((2-(2-((6-(bis(carboxylatomethyl)amino)-2-(5-carboxylatooxazol-2-yl)benzofuran-5-yl)oxy)ethoxy)-4-methylphenyl)azanediy)diacetate, pentasodium salt

**MF:** C<sub>29</sub>H<sub>22</sub>N<sub>3</sub>O<sub>14</sub> • 5Na

**FW:** 751.5

**Purity:** ≥90%

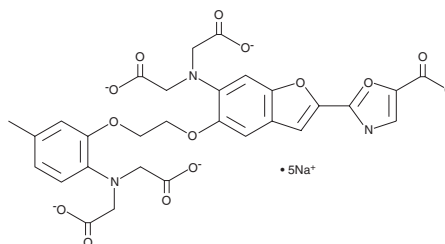
**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥2 years

**Special Conditions:** Protect from light

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.



### Laboratory Procedures

Fura-2 (sodium salt) is supplied as a solid. It is soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Fura-2 is a ratiometric fluorescent calcium indicator that can be used to detect calcium in cells.<sup>1</sup> It is a pentacarboxylate that displays excitation maxima of 340 and 380 nm at high and low calcium concentrations, respectively, when the emission is fixed at 510 nm, enabling determination of ratiometric measurements of calcium influx in live cells.

### Reference

1. Kong, S.K. and Lee, C.Y. The use of fura 2 for measurement of free calcium concentration. *Biochem. Mol. Biol. Ed.* **23(2)**, 97-98 (1995).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/06/2019

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM