

PRODUCT INFORMATION



Fura-2 Leakage Resistant (potassium salt)

Item No. 34994

Formal Name: 2,2'-((2-(2-((6-(bis(carboxylatomethyl)amino)-2-(5-carboxylatooxazol-2-yl)benzofuran-5-yl)oxy)ethoxy)-4-(2-carboxylatoethyl)phenyl)azanediyl) diacetate, pentapotassium salt

Synonyms: Fura-2 LR, Fura-PE3

MF: $C_{31}H_{23}K_6N_3O_{16} \bullet 5K$

FW: 1,123.6

Purity: $\geq 90\%$

Ex./Em. Max: 340 and 380/505 nm

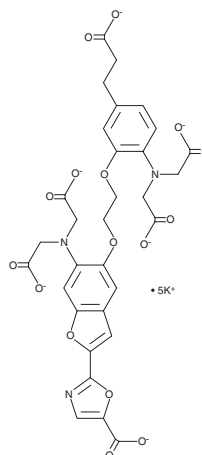
Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

Special Conditions: Protect from light and moisture

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



Laboratory Procedures

Fura-2 leakage resistant (potassium salt) is supplied as a solid. A stock solution may be made by dissolving the fura-2 leakage resistant (potassium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Fura-2 leakage resistant is a cell-impermeable, ratiometric fluorescent calcium indicator. It binds to calcium ($K_d = 145$ nM) and displays excitation maxima of 340 and 380 nm when calcium-bound or calcium-free, respectively, when the emission is fixed at 505 nm. Fura-2 leakage resistant can be used in lipid membrane-free systems or liposomes or can be introduced into cells by electroporation, microinjection, or other methods. Fura-2 leakage resistant is also available in a cell-permeable form (Item No. 34993).

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, 06/05/2024

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