

PRODUCT INFORMATION



ICR-1 (potassium salt)

Item No. 34986

Synonym: ION Calcium Red-1
Purity: $\geq 90\%$
UV/Vis.: λ_{max} : 580 nm
Ex./Em. Max: 580/660 nm
Supplied as: A solid
Storage: -20°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.

Description

ICR-1 is a cell-impermeable fluorescent calcium indicator. It binds to calcium ($K_d = 480$ nM) and displays excitation/emission maxima of 580/660 nm, respectively. ICR-1 can be used in lipid membrane-free systems or liposomes or can be introduced into cells by electroporation, microinjection, or other methods. ICR-1 is also available in a cell-permeable form (Item No. 34985).

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, 12/08/2022

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