

PRODUCT INFORMATION



Fluo-2 (potassium salt)

Item No. 35764

Formal Name: N-[2-[2-[2-[bis(carboxymethyl)amino]-5-(6-hydroxy-3-oxo-3H-xanthen-9-yl)phenoxy]ethoxy]-4-methylphenyl]-N-(carboxymethyl)glycine, pentapotassium salt

Synonyms: Fluo-2 HA, Fluo-2 High Affinity

MF: C₃₆H₂₇N₂O₁₃ • 5K

FW: 891.1

Purity: ≥95%

UV/Vis.: λ_{max}: 254 nm

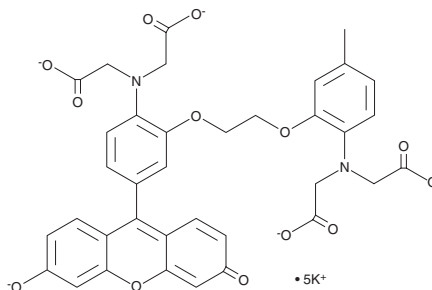
Ex./Em. Max: 490/515 nm

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

Fluo-2 is a cell-impermeable fluorescent calcium indicator. It binds to calcium ($K_d = 290$ nM) and displays excitation/emission maxima of 490/515 nm, respectively. Fluo-2 is also available in a cell-permeable form (Item No. 35758).

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, 03/29/2024

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