

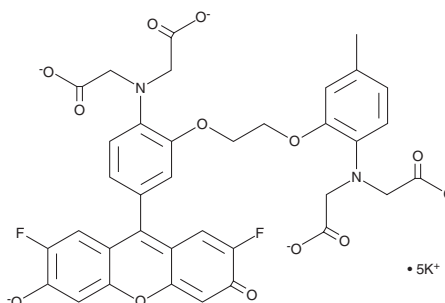
PRODUCT INFORMATION



Fluo-4 (potassium salt)

Item No. 35765

CAS Registry No.: 1199351-32-0
Formal Name: N-[2-[2-[2-[bis(carboxymethyl)amino]-5-(2,7-difluoro-6-hydroxy-3-oxo-3H-xanthen-9-yl)phenoxy]ethoxy]-4-methylphenyl]-N-(carboxymethyl)glycine, pentapotassium salt
MF: C₃₆H₂₅F₂N₂O₁₃ • 5K
FW: 927.1
Purity: ≥90%
Ex./Em. Max: 490/515 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.

Laboratory Procedures

Fluo-4 (potassium salt) is supplied as a solid. A stock solution may be made by dissolving the fluo-4 (potassium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Fluo-4 is a cell-impermeable fluorescent calcium indicator. It binds to calcium ($K_d = 355$ nM) and displays excitation/emission maxima of 490/515 nm, respectively.

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