

PRODUCT INFORMATION



Fluo-3FF (potassium salt)

Item No. 20406

CAS Registry No.: 348079-14-1
Formal Name: N-[2-[2-[2-[bis(carboxymethyl)amino]-5-(2,7-dichloro-6-hydroxy-3-oxo-3H-xanthen-9-yl)phenoxy]ethoxy]-3,4-difluorophenyl]-N-(carboxymethyl)-glycine, pentapotassium salt

MF: C₃₅H₂₁Cl₂F₂N₂O₁₃ • 5K

FW: 981.9

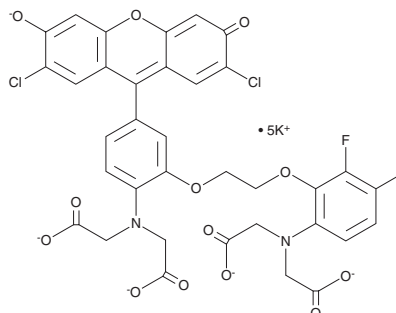
Purity: ≥90%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Special Conditions: Protect from light



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

Laboratory Procedures

Fluo-3FF (potassium salt) is supplied as a solid. A stock solution may be made by dissolving the fluo-3FF (potassium salt) in the solvent of choice. Fluo-3FF (potassium salt) is soluble in water.

Description

Fluo-3FF (potassium salt) is a fluorescent calcium indicator. It is a di-fluorinated analog of Fluo-3 (Item Nos. 20402 | 20403 | 20404) that displays about a 100-fold lower affinity than fluo-3 for calcium ($K_{dS} = 42$ and $0.4 \mu\text{M}$, respectively).¹ Low affinity calcium indicators are particularly useful for studying compartments with high concentrations of calcium, such as endoplasmic reticulum, where high affinity dyes will be insensitive to luminal fluctuations. Fluo-3FF has excitation/emission maxima of approximately 506 and 526 nm, respectively.

Reference

1. David, G., Barrett, J. N. and Barrett, E. F. Stimulation-induced changes in Ca^{2+} in lizard motor nerve terminals. *J. Physiol.* **504**(Pt 1), 83-96 (1997).

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, 11/03/2022

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