

PRODUCT INFORMATION



Rhod-FF (potassium salt)

Item No. 20443

Formal Name: 2,2'-((2-(2-(6-(bis(carboxylatomethyl)amino)-2,3-difluorophenoxy)ethoxy)-4-(6-(dimethylamino)-3-(dimethyliminio)-3H-xanthen-9-yl)phenyl)azanediyl) diacetate, tripotassium salt

MF: C₃₉H₃₅F₂N₄O₁₁ • 3K

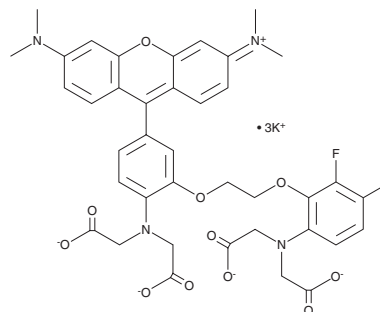
FW: 891.0

Purity: ≥90%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

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

Description

Rhod-FF is a difluorinated analog of the cell-impermeant calcium indicator rhod-2. It has a very low affinity for calcium ($K_d = 320 \mu\text{M}$).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. Rhod-FF has excitation/emission maxima of approximately 552 and 580 nm, respectively.

Reference

1. Paredes, R. M., Etzler, J. C., Watts, L. T., *et al.* Chemical calcium indicators. *Methods* **46(3)**, 143-151 (2008).

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, 10/18/2022

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