

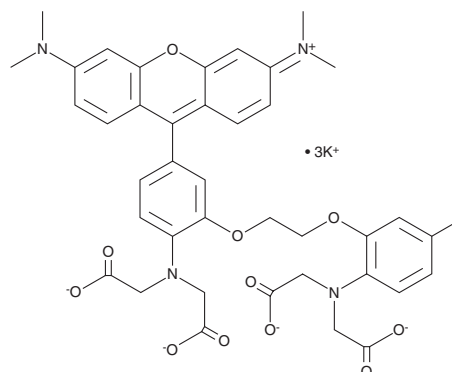
PRODUCT INFORMATION



Rhod-2 (potassium salt)

Item No. 20777

CAS Registry No.: 663625-18-1
Formal Name: 2,2'-((2-(2-(2-(bis(carboxylato methyl)amino)-5-(6-(dimethylamino)-3-(dimethyliminio)-3H-xanthen-9-yl) phenoxy)ethoxy)-4-methylphenyl) azanediyl)diacetate, tripotassium salt
MF: C₄₀H₃₉N₄O₁₁ • 3K
FW: 869.1
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Avoid light



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

Laboratory Procedures

Rhod-2 (potassium salt) is supplied as a solid. A stock solution may be made by dissolving the Rhod-2 (potassium salt) in the solvent of choice, which should be purged with an inert gas. Rhod-2 (potassium salt) is soluble in DMSO. It is also soluble in water.

Description

Rhod-2 (potassium salt) is a water-soluble, red fluorescent calcium indicator. It exhibits a significant shift in fluorescence intensity upon calcium binding (ex max = 549 nm; calcium-free v. ex/em max = 552/581 nm; calcium-bound).^{1,2} Unlike the UV-excitable indicators fura-2 (Item Nos. 19531 | 20414) and indo-1 (Item Nos. 20419 | 20418), there is no accompanying spectral shift.

References

1. Paredes, R.M., Etzler, J.C., Watts, L.T., *et al.* Chemical calcium indicators. *Methods: A Companion to Methods in Enzymology* **46(3)**, 143-151 (2008).
2. Minta, A., Kao, J.P., Tsien, R.Y. Fluorescent indicators for cytosolic calcium based on rhodamine and fluorescein chromophores. *Journal of Biological Chemistry* **264(14)**, 8171-8178 (1989).

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, 10/21/2022

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