

PRODUCT INFORMATION



Rhod-2 (sodium salt)

Item No. 20778

Formal Name: 2,2'-((2-(2-(2-(bis(carboxylatomethyl)amino)-5-(6-(dimethylamino)-3-(dimethyliminio)-3H-xanthen-9-yl)phenoxy)ethoxy)-4-methylphenyl)azanediyl)diacetate, trisodium salt

MF: C₄₀H₃₉N₄O₁₁ • 3Na

FW: 820.7

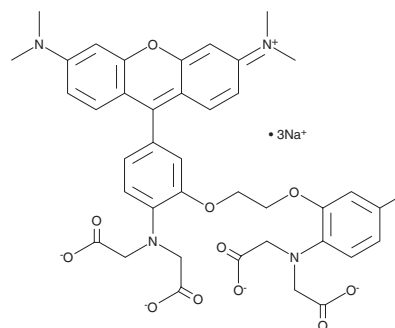
Purity: ≥90%

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 (sodium salt) is supplied as a solid. It is moderately soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Rhod-2 (sodium 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, 11/12/2022

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