## Rhod-2, Tripotassium salt

Membrane-impermeant calcium indictor that can be loaded into cells via microinjection or scrape loading.



## Product attributes

Cell permeability	Membrane impermeant
,	
Indicator type	Non-ratiometric
Excitation/Emission	556/576 nm

Call us: 800-304-5357 Email: btinfo@biotium.com

## **Product Description**

Membrane-impermeant calcium indictor that can be loaded into cells via microinjection or scrape loading.

- $\lambda_{Ex}/\lambda_{Em}$  (low or high [Ca<sup>2+</sup>])= 556/576 nm
- $\epsilon$  (556 nm) = 80,000 M<sup>-1</sup>cm<sup>-1</sup>.
- Orange red solid soluble in DMSO and water (pH >6)
- Store at 4°C and protect from light, especially in solution
- C40H39K3N4O11
- MW: 869

BAPTA-based ion indicators like Rhod-2 have been shown to be fixable in situ by EDC/EDAC (cat# 59002). The fixation of indicator dyes is useful for downstream immunofluorescence and IHC studies (Cell Calcium 1997, 21(3), 175).

As the indicator does not covalently bind to cellular components, it may be actively effluxed from the cell by organic anion transporters. The rate of efflux increases with temperature, and may vary between cell types, resulting in variable retention times of a few minutes to hours. Experiments using indicators in cells usually are performed within one or two hours of loading, but it may be possible to re-load cells with indicator if needed. The organic anion transporter inhibitor Probenecid (#50027) can be used to slow the rate of indicator efflux from cells.

## References

- 1. Biophys J, 59, 12 (1991), <u>DOI: 10.1016/S0006-3495(91)82193-2</u>
  2. Science 250, 977 (1990), <u>DOI: 10.1126/science.2237441</u>
  3. J Biol Chem 264, 8171 (1989), <u>DOI: 10.1016/S0021-9258(18)83165-9</u>
- 4. Methods Cell Biol, 99, 113, (2021), DOI: 10.1016/B978-0-12-374841-6.00005-0

This datasheet was generated on November 9, 2025 at 08:24:19 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/rhod-2-tripotassium-salt/