RH795

The styryl dye RH795 is a fast-responding potentiometric probe that is primarily used for functional imaging of neurons. The dye is spectrally similar to RH414 (61016), but exhibits different physiological effects during staining. For example, RH414 causes arterial constriction during cortex staining, while RH795 does not (see references 1,2).



Product attributes

CAS number	172807-13-5
For live or fixed cells	For live/intact cells
Potential dependence	Fast-response membrane potential dye
Colors	Far-red
Excitation/Emission	530/712 nm (see product description)

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

Product Description

The styryl dye RH795 is a fast-responding potentiometric probe that is primarily used for functional imaging of neurons. The dye is spectrally similar to RH414 (#61016), but exhibits different physiological effects during staining. For example, RH414 causes arterial constriction during cortex staining, while RH795 does not (see references 1,2). The excitation/emission data below are for the dye in methanol. In cell membranes, the spectra of styryl dyes are typically blue-shifted by as much as 20 nm for absorption or excitation and 80 nm for emission.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 530/712 nm
- Orange red solid soluble in H2O
- Store at -20°C and protect from light
- C₂₆H₃₉Br₂N₃O₂
- MW: 585.42
- [172807-13-5]

References

- 1. Nature 324, 361 (1986).
- 2. J Neurosci 14, 2545 (1994).
- 3. PLoS ONE 8(10), e75678 (2013).

This datasheet was generated on November 8, 2025 at 11:56:45 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/rh795/