

RH237

The styryl dye RH237 is a fast-responding potentiometric probe that is primarily used for functional imaging of neurons.



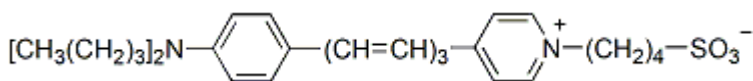
Product attributes

For live or fixed cells	For live/intact cells
Potential dependence	Fast-response membrane potential dye
Colors	Far-red
Excitation/Emission	528/782 nm (see product description)

Product Description

The styryl dye RH237 is a fast-responding potentiometric probe that is primarily used for functional imaging of neurons. Excitation/emission data (below) is 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) = 528/782 nm
- Dark solid soluble in DMSO
- Store at -20 °C and protect from light
- $C_{29}H_{40}N_2O_3S$
- MW: 496.71



References

1. Neuroscience 31, 613 (1989).

This datasheet was generated on January 11, 2026 at 08:20:39 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/rh237/>