

RH237

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



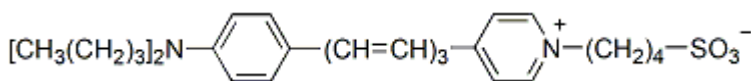
Product attributes

| | |
|-------------------------|--------------------------------------|
| 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 June 12, 2026 at 10:05:09 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/rh237/>