Glowing products for science

RH414

The styryl dye RH414 is a fast-responding potentiometric probe that is primarily used for functional imaging of neurons.



Product attributes

CAS number	161433-30-3
For live or fixed cells	For live/intact cells
Potential dependence	Fast-response membrane potential dye
Colors	Far-red
Excitation/Emission	532/716 nm (see product description)

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

Product Description

The styryl dye RH414 is a fast-responding potentiometric probe that is primarily used for functional imaging of neurons. Excitation/emission data 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) = 532/716 nm
- Orange red soluble in H2O or DMSO
- Store at -20°C and protect from light
- C₂₈H₄₃Br₂N₃
- MW: 581.48
- [161433-30-3]

$$2Br^{-}$$
 $(CH_{3}CH_{2})_{2}N$
 $(CH=CH)_{2}$
 $N^{+}(CH_{2})_{3}N^{+}(CH_{2}CH_{3})_{3}$

References

1. Neuroscience 31, 613 (1989).

This datasheet was generated on November 2, 2025 at 05:08:38 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/rh414/