

## RH414

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



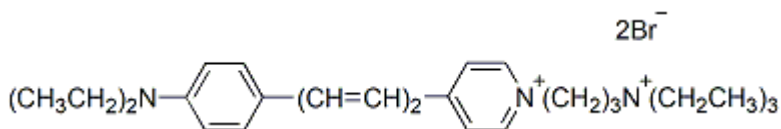
### Product attributes

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

## 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 H<sub>2</sub>O or DMSO
- Store at -20 °C and protect from light
- C<sub>28</sub>H<sub>43</sub>Br<sub>2</sub>N<sub>3</sub>
- MW: 581.48
- [161433-30-3]



## References

1. Neuroscience 31, 613 (1989).

This datasheet was generated on June 22, 2026 at 11:18:22 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/rh414/>