Octadecyl Rhodamine B Chloride (R18)

The fluorescence of octadecyl rhodamine B in membranes is quenched at high dye concentration but is released at dilution. This property makes the dye useful for membrane fusion assays.



Product attributes

CAS number	65603-19-2
Probe cellular localization	Membrane/cell surface, Membrane/vesicular
Colors	Red
Excitation/Emission	556/578 nm

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

Product Description

The fluorescence of octadecyl rhodamine B in membranes is quenched at high dye concentration but is released at dilution. This property makes the dye useful for membrane fusion assays.

- λEx/λEm (MeOH) = 556/578 nm
- Dark solid soluble in EtOH or DMSO
- Store at -20°C and protect from light
- C46H67CIN2O3
- MW: 731.50
- [65603-19-2]

This datasheet was generated on November 4, 2025 at 09:54:22 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/octadecyl-rhodamine-b-chloride-r18/