

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.



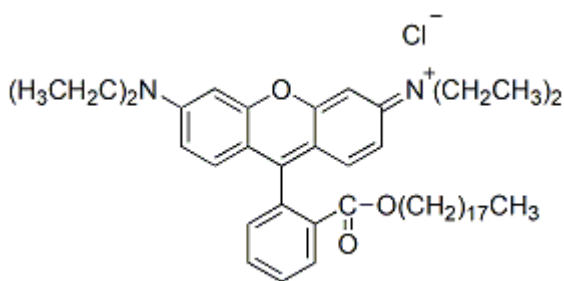
Product attributes

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

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.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 556/578 nm
- Dark solid soluble in EtOH or DMSO
- Store at -20 °C and protect from light
- $C_{46}H_{67}ClN_2O_3$
- MW: 731.50
- [65603-19-2]



This datasheet was generated on June 4, 2026 at 06:46:09 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/octadecyl-rhodamine-b-chloride-r18/>