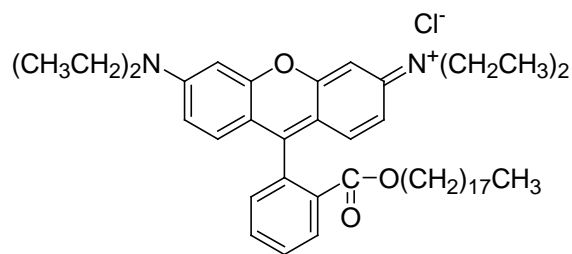


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Octadecyl rhodamine B chloride (R18)**CATALOG #:** 60033**MOLECULAR INFORMATION:** C<sub>46</sub>H<sub>67</sub>C<sub>1</sub>N<sub>2</sub>O<sub>3</sub>  
Mwt: 731.5**PROPERTIES:**

<b>Color &amp; Form</b>	Red solid
<b>Purity</b>	≥ 95% by TLC
<b>Solubility</b>	Soluble in DMSO or EtOH
$\lambda_{ex} \backslash \lambda_{em}$	556 nm/578 nm (MeOH)
<b>Extinction Coefficient</b>	100,000 (MeOH)

**STORAGE AND HANDLING:** Store at -20 °C and protect from light especially when in solution.**APPLICATION:** 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.**TOXICITY:** Unknown.**FIRST AID:** Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.**Disclaimer:** Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.