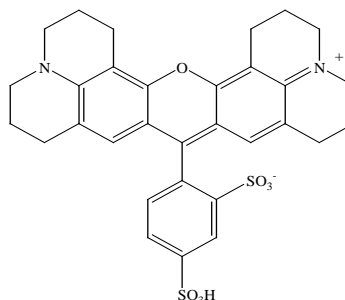


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Sulforhodamine 101**CATALOG #:** 80101**MOLECULAR INFORMATION:** C<sub>31</sub>H<sub>30</sub>N<sub>2</sub>O<sub>7</sub>S<sub>2</sub>  
Mwt: 607  
[60311-02-6]**PROPERTIES:**

<b>Color &amp; Form</b>	Red solid
<b>Purity</b>	98% by HPLC
<b>Solubility</b>	Soluble in water
<b>Absorption/Emission</b>	576/605 nm
<b>Extinction Coefficient</b>	>80,000

**STORAGE AND HANDLING:** Store at room temperature and protect from light.**APPLICATION:** Sulforhodamine 101 has a longer absorption wavelength than does sulforhodamine B. Like sulforhodamine B and G, in addition to their potential use in cancer drug screening, these fluorescent dyes have been primarily used as polar tracers for the studies of neuronal cell morphology and cell-cell communications<sup>1</sup>. See **80100** for more information.

Ref: 1: a) Meth. Enzymol. 221, 234(1993); b) J. Biol. Chem. 267, 18424(1992); c) Plant Cell Environ.17, 257(1994)

**TOXICITY:** Unknown.

<b>FIRST AID:</b>	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.
-------------------	---

<b>Disclaimer:</b> 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.
---