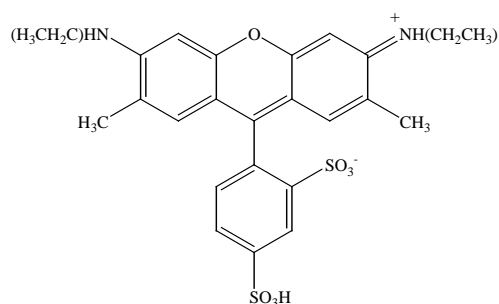


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** **Sulforhodamine G****CATALOG #:** **80102****MOLECULAR INFORMATION:** C<sub>25</sub>H<sub>25</sub>N<sub>2</sub>NaO<sub>7</sub>S<sub>2</sub>  
MWt: 553  
[5873-16-5]**PROPERTIES:**

<b>Color &amp; Form</b>	Orange red solid
<b>Purity</b>	≥ 95% by TLC
<b>Solubility</b>	Soluble in water
<b>Absorption/Emission</b>	532/548 nm
<b>Extinction Coefficient</b>	

**STORAGE AND HANDLING:**

Store at room temperature. Protect material from light, especially in solution.

**APPLICATION:**

Sulforhodamine G has a shorter absorption wavelength. 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.

Biotium also offers **sulforhodamine 101(80101)** and **sulforhodamine B(80100)**. 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.

**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> <i>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.</i>
--