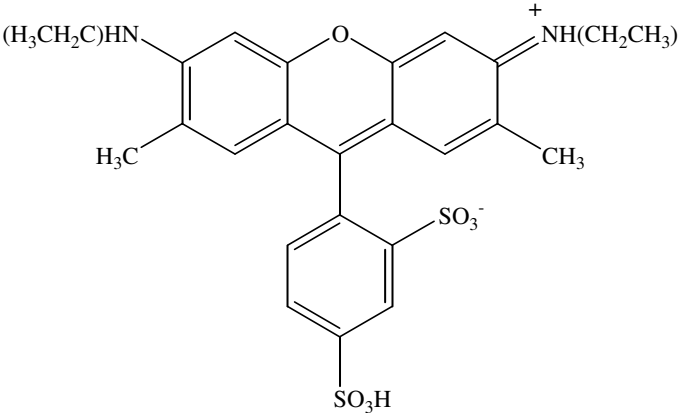


Instruction Manual

Catalog Number	PK-CA707-80102
Description	Similar to other sulforhodamines, the highly water-soluble sulforhodamine G can also be used as a polar tracer for the studies of cell morphology and neuronal cell-cell communications.
Quantity	5 g
Excitation / Emission Maxima	$\lambda_{\text{exc}} \setminus \lambda_{\text{em}}$: 532/548 nm
Molecular Structure	
Molecular Weight / Molecular Formula	553 Da; C ₂₅ H ₂₅ N ₂ NaO ₇ S ₂
Purity	>95% (as determined by TLC)
Appearance / Formulation / Solubility	Orange-red solid; soluble in water.
Storage & Stability	Store at room temperature. Protect material from light, especially when in solution.
Applications	See Description
References	<ol style="list-style-type: none"> 1) Meth. Enzymol. 221, 234(1993) 2) J. Biol. Chem. 267, 18424(1992) 3) Plant Cell Environ. 17, 257(1994)
Caution	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.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.