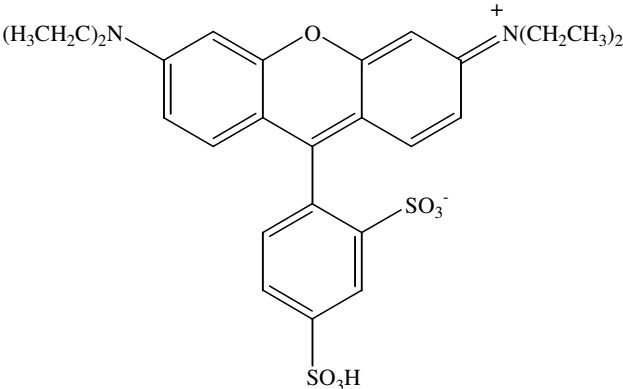


Instruction Manual

Catalog Number	PK-CA707-80100
Description	Similar to other sulforhodamines, the highly water-soluble sulforhodamine B can also be used as a polar tracer for the studies of cell morphology and neuronal cell-cell communications. In addition, sulforhodamine B has been shown to be a superior protein stain for use in the quantification of cellular proteins of cultured cells. The intensely colored, water soluble dye is particularly useful for in vitro cell-based screening of anticancer drugs. The dye is believed to bind to basic amino acids of cellular proteins. Thus, colorimetric measurement of the bound dye provides an estimate of the total protein mass that is related to the cell number. The assay method is simple and reproducible, and the end-point measurement is not time-critical, a significant advantage over assays using tetrazolium derivatives. The method is especially applicable to large scale screenings of anticancer drug candidates from natural extracts or synthetic chemicals.
Quantity	5 g
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em} = 565/586 \text{ nm}$
Molecular Structure	
Molecular Weight / Molecular Formula	559 Da; C ₂₇ H ₃₀ N ₂ O ₇ S ₂
Purity	>95% (as determined by TLC)
Appearance / Formulation / Solubility	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) 4) J. Natl. Cancer Inst. 82, 1107(1990) 5) J. Natl. Cancer Inst. 82, 1087(1990) 6) J. Immunol. Meth. 208, 151(1997)
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.