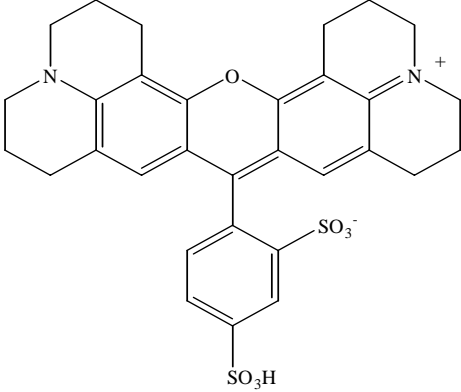


Instruction Manual

Catalog Number	PK-CA707-80101
Description	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. See PK-CA707-80100 for more information.
Quantity	500 mg
Excitation / Emission Maxima	λ_{ex} \ λ_{em} : 586/605 nm; Extinction coefficient = 108,000
Molecular Structure	
Molecular Weight / Molecular Formula	607 Da; C ₃₁ H ₃₀ N ₂ O ₇ S ₂
Purity	>98% (as determined by HPLC)
Appearance / Formulation / Solubility	Red solid; soluble in water.
Storage & Stability	Store at room temperature. Protect material from light.
Applications	See Description
References	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.