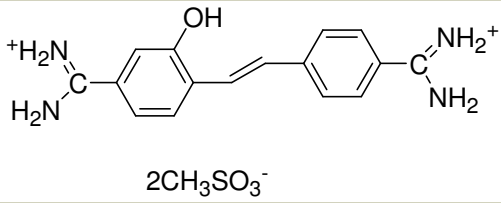


Instruction Manual

Catalog Number	PK-CA707-80023
Description	Hydroxystilbamidine has been used extensively as a retrograde tracer for neurons and also a histochemical stain. The 4% solution in deionized water is a more convenient, ready-to-use form. If desired, the solution can be diluted to a lower concentration. For most of the applications, the typically dye concentration is 2% to 4%. Hydroxystilbamidine has very broad spectra at physiological pH with absorption from 350nm to 400nm and emission from 530 to 620 nm. A wide band UV excitation filter is recommended for optimal excitation.
Quantity	200 µl
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em} = 361/536$ nm. λ_{abs} : 290 ~ 370nm (broad absorption in water). When bound to membranes at pH 7.4, excitation band: 350-395 nm (broad peak); emission band: 530-600 nm (broad peak).
Molecular Structure	 <p style="text-align: center;">$2\text{CH}_3\text{SO}_3^-$</p>
Molecular Weight / Molecular Formula	473 Da; $\text{C}_{18}\text{H}_{24}\text{N}_4\text{O}_7\text{S}_2$
Purity	>95% (as determined by HPLC)
Appearance / Formulation / Solubility	Yellow liquid; soluble in water.
Storage & Stability	Store desiccated at 4°C. Protect from light.
Applications	See Description
References	1) Brain Res. 377, 147(1986) 2) Brain Res. 475, 259(1988) 3) J. Neurosci. methods, 23, 1(1988); 4) Brain Res. 553, 135(1991) 5) J. Neurosci. methods, 45, 87(1992)
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.

*FluoroGold is a trademark from Fluorochrome, LLC

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