

RH 795

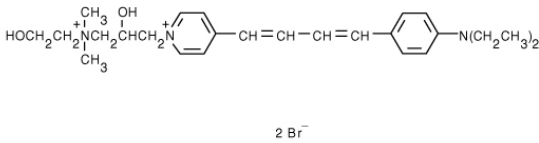
Catalog Number	Packaging Size
C267	1 mg

Storage upon receipt: -20°C, protected from light

Introduction

RH 795 is a dialkylaminophenylpolyenylpyridinium dye that is principally used for functional imaging of neurons. RH 795 produced negligible side effects when tested in vitro using hippocampal slices and in vivo using single-unit recordings in cat and monkey visual cortices.

Specifications

Label:	RH 795	
Ex/Em:	486/689 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula:	C ₂₆ H ₃₉ Br ₂ N ₃ O ₂	
Molecular Weight:	585.42	
CAS Number:	172807-13-5	
Storage Conditions:	-20°C, protected from light	
Shipping Condition:	Room Temperature	

Applications

Membrane potential indicator

References:

[A new neurophysiological/neuropathological ex vivo model localizes the origin of glioma-associated epileptogenesis in the invasion area.](#)

Senner V, Köhling R, Püttmann-Cyrus S, Straub H, Paulus W, Speckmann EJ
Acta Neuropathol (Berl) (2004) 107:1-7

[Novel naphthylstyryl-pyridium potentiometric dyes offer advantages for neural network analysis.](#)

Obaid AL, Loew LM, Wuskell JP, Salzberg BM
J Neurosci Methods (2004) 134:179-190

[Optical current source density analysis in hippocampal organotypic culture shows that spreading depression occurs with uniquely reversing currents.](#)

Kunkler PE, Hulse RE, Schmitt MW, Nicholson C, Kraig RP
J Neurosci (2005) 25:3952-3961