

Product Information Sheet

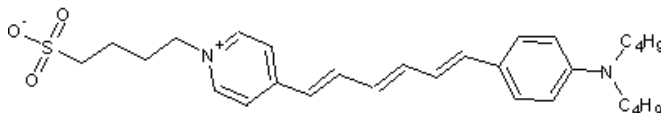
Ordering Information

Product Number: 21480
Product Name: RH 237
Unit Size: 5 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 496.7

Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 528 nm
Emission Wavelength: 782 nm

Application Notes

A neuron-tracing dye. It is used for monitoring membrane potential, synaptic activity and ion channel activity of neurons.