

## RH 237 [N-(4-Sulfobutyl)-4-(6-(4-(dibutylami no)phenyl)hexatrienyl)pyridinium, inner salt]

Catalog number: 21480 Unit size: 5 mg

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties	
Appearance	Dark brown solid
Molecular Weight	496.7
Soluble In	DMSO
Chemical Structure	

## Applications

RH 237, also called N-(4-sulfobutyl)-4-(6-(4-(dibutylamino)phenyl)hexatrienyl)pyridinium, is a neuron-tracing dye. It is used for monitoring membrane potential, synaptic activity and ion channel activity of neurons.