

**RH 155 \*CAS#: 254451-42-8\***

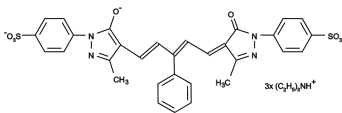
Catalog number: 21493

Unit size: 25 mg

## Product Details

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

## Chemical Properties

Appearance	Solid
Molecular Weight	950.26
Soluble In	Water
Chemical Structure	

## Applications

RH 155 is a neuron-tracing dye. It is used for monitoring membrane potential, synaptic activity and ion channel activity of neurons.