

RH 414 [N-(3-Triethylammoniumpropyl)-4-(4-(4-(diethylamino)phenyl)butadienyl)pyridi nium dibromide] *CAS#: 161433-30-3*

Catalog number: 21485

Unit size: 5 mg

Product Details

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Dark blue solid

Molecular Weight 581.47

Soluble In Water

Chemical Structure

Applications

RH 414 [N-(3-Triethylammoniumpropyl)-4-(4-(diethylamino)phenyl)butadienyl)pyridinium dibromide] is a neuron-tracing dye. It is used for monitoring membrane potential, synaptic activity and ion channel activity of neurons.