

Product Information

Dihydrorhodamine 123

Catalog Number	Unit Size
C261	10 mg

Storage upon receipt:

- -20°C
- Protected from light

Product Description

Dihydrorhodamine 123 is an uncharged and nonfluorescent reactive oxygen species (ROS) indicator that can passively diffuse across membranes where it is oxidized to cationic rhodamine 123 which localizes in the mitochondria and exhibits green fluorescence.

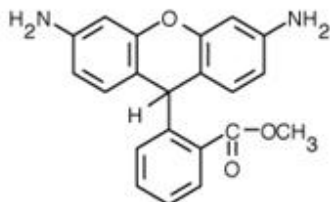
Specifications

Chemical Name: Dihydrorhodamine 123

Molecular Formula: C₂₁H₁₈N₂O₃

Molecular Weight: 346.38

CAS Number: 109244-58-8



Application

Reactive oxygen species (ROS) indicator.