

DHR 6G [Dihydrorhodamine 6G]

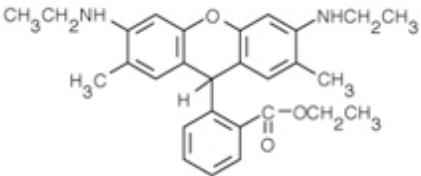
Catalog Number	Packaging Size
C262	25 mg

Storage upon receipt: -20°C, protected from light

Introduction

Dihydrorhodamine 6G is the reduced form of rhodamine 6G, which is used as fluorescent mitochondrial dye. It is nonfluorescent, but it readily enters most of the cells and is oxidized by oxidative species or by cellular redox systems to the fluorescent rhodamine 6G that accumulates in mitochondrial membranes. Dihydrorhodamine 6G is useful for detecting reactive oxygen species (ROS) including superoxide.

Specifications

Label:	Rhodamine 6G	
Ex/Em:	528/551 nm	
Detection Method:	Fluorescent	
Molecular Formula:	C ₂₈ H ₃₂ N ₂ O ₃	
Molecular Weight:	444.57	
CAS Number:	217176-83-5	
Storage Conditions:	-20°C, protected from light	
Shipping Condition:	Room Temperature	

Applications

Probe for ROS

References:

1. [Fluorescent approach to quantitation of reactive oxygen species in mainstream cigarette smoke.](#)
Ou B, Huang D
Anal Chem (2006) 78:3097-3103
2. [Uptake of fluorescent dyes associated with the functional expression of the cystic fibrosis transmembrane conductance regulator in epithelial cells.](#)
Wersto RP, Rosenthal ER, Crystal RG, Spring KR
Proc Natl Acad Sci U S A (1996) 93:1167-1172
3. [Kindlin-3 is required for beta2 integrin-mediated leukocyte adhesion to endothelial cells.](#)
Moser M, Bauer M, Schmid S, Ruppert R, Schmidt S, Sixt M, Wang HV, Sperandio M, Fässler R,
Nat Med (2009) 15:300-305