

Resorufin β -D-galactopyranoside

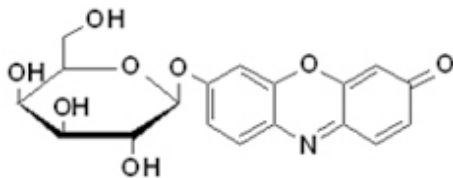
Catalog Number	Packaging Size
C280	10 mg

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

Introduction

The fluorogenic β -galactosidase substrate **resorufin β -D-galactopyranoside** yields the hydrolysis product, resorufin, with absorption/emission maxima at 571/585 nm.

Specifications

Label:	Resorufin	
Ex/Em:	571/585 nm	
Detection Method:	Fluorescent	
Molecular Formula:	C ₁₈ H ₁₇ NO ₈	
Molecular Weight:	375.33	
CAS Number:	95079-19-9	
Storage Conditions:	-20°C, protected from light	
Shipping Condition:	Room Temperature	

Applications

Galactosidase Substrate

References:

- Flow Fluorimetric Investigations on Reaction Kinetics of a Growing Analytical Bioreactor.
Willems H, Sernetz M
Anal Chim Acta (1988) 213:245-245
- Immobilized Enzyme Kinetics Analyzed by Flow-Through Microfluorimetry.
Resorufin-B-D-galactopyranoside as a New Fluorogenic Substrate for Galactosidase.
Hofmann J, Sernetz M
Anal Chim Acta (1984) 163:67-67
- [Subnanoliter enzymatic assays on microarrays.](#)
Angenendt P, Lehrach H, Kreutzberger J, Glökler J
Proteomics (2005) 5:420-425