

CUG (3-Carboxyumbelliferyl β-D-galactopyranoside)

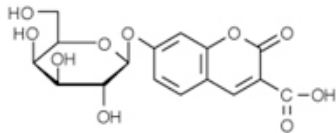
Catalog Number	Packaging Size
C277	10 mg

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

Introduction

CUG (3-Carboxyumbelliferyl β-D-galactopyranoside) is a β-galactosidase substrate that produces the blue-fluorescent hydrolysis product, 7-hydroxycoumarin-3-carboxylic acid.

Specifications

Label:	7-hydroxycoumarin-3-carboxylic acid	
Ex/Em:	386/448 nm	
Detection Method:	Fluorescent	
Molecular Formula:	C ₁₆ H ₁₆ O ₁₀	
Molecular Weight:	368.30	
CAS Number:	64664-99-9	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Galactosidase Substrate