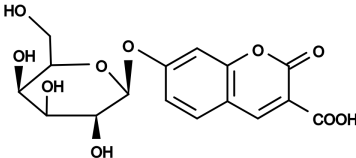


CUG [3-Carboxyumbelliferyl beta-D-galactopyranoside] *CAS 64664-99-9*Catalog number: 14003
Unit size: 10 mg**Product Details**

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	368.29
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	352 nm
Emission Wavelength	407 nm

Applications

CUG is a fluorogenic galactosidase substrate with carboxy reactive handle that can be used for conjugation. The organic and inorganic esters and glycosides of 7-hydroxycoumarin are among the most commonly used fluorogenic substrates. These substrates yield water-soluble blue fluorescent coumarin fluorophores upon interaction with enzymes.