

Product Information Sheet

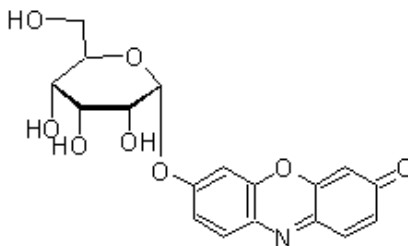
Ordering Information

Product Number: 14021
Product Name: Resorufin alpha-D-galactopyranoside
Unit Size: 5 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 375.33

Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 571 nm
Emission Wavelength: 585 nm

Application Notes

Resorufin alpha-D-galactopyranoside is a sensitive fluorogenic substrate that generates a red fluorescent product (resorufin) upon interaction with alpha-galactosidase. It is used for measuring alpha-galactosidase activities and high throughput screening of alpha-galactosidase inhibitors. Alpha-galactosidase is a glycoside hydrolase enzyme that hydrolyses the terminal alpha-galactosyl moieties from glycolipids and glycoproteins. This enzyme is a homodimeric glycoprotein that hydrolyses the terminal alpha-galactosyl moieties from glycolipids and glycoproteins. It predominantly hydrolyzes ceramide trihexoside, and it can catalyze the hydrolysis of melibiose into galactose and glucose. A variety of mutations in this gene affect the synthesis, processing, and stability of this enzyme, which causes Fabry's disease, a rare lysosomal storage disorder and sphingolipidosis that results from a failure to catabolize alpha-D-galactosyl glycolipid moieties.