

Resorufin alpha-D-galactopyranoside

Catalog number: 14021

Unit size: 5 mg

Product Details

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Orange solid

Molecular Weight 375.33

Soluble In DMSO

Chemical Structure

Spectral Properties

Excitation Wavelength 571 nm

Emission Wavelength 584 nm

Applications

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.