

Product Information Sheet

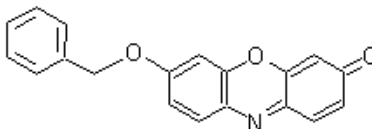
Ordering Information

Product Number:	15020
Product Name:	Resorufin benzyl ether
Unit Size:	10 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 303.31

Chemical Structure:



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

Application Notes

This fluorogenic resorufin ether is widely used for monitoring P450 activities in cell extracts and solutions. Cytochrome P450 oxidase (CYP450) is a large number of evolutionary related oxidative enzymes important in animal, plant, and bacterial physiology. Most cytochromes P450 (CYPs) have about 500 amino acids and a heme group at the active site.