

## Product Information Sheet

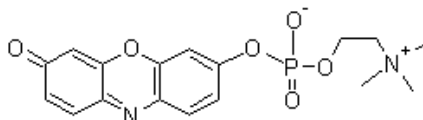
### Ordering Information

Product Number:	11640
Product Name:	Resorufin Phosphocholine
Unit Size:	1 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

### Chemical and Spectral Properties

Molecular Weight: 378.32

Chemical Structure:



Soluble in:	Water
Excitation Wavelength:	575 nm
Emission Wavelength:	590 nm

### Application Notes

Glycerophosphocholine phosphodiesterase release phosphocholine from resorufin phosphocholine to give the highly fluorescent resorufin that can be readily monitored by either fluorescence or absorption. This fluorogenic phosphodiesterase substrate can be conveniently used for screening the inhibitors of glycerophosphocholine phosphodiesterase. It can also be used for monitoring the activities of phospholipase D in combination with a phosphatase. Phospholipase D hydrolyzes resorufin phosphocholine to resorufin phosphate, which in the presence of an acid phosphatase generates resorufin. Thus, resorufin phosphocholine can also be used for monitoring phospholipase D.