

Resorufin Phosphocholine

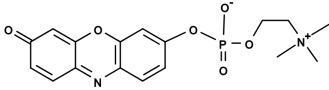
Catalog number: 11640

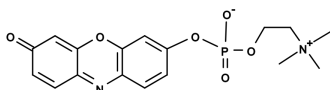
Unit size: 1 mg

Product Details

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

Chemical Properties

Appearance	Yellow solid
Molecular Weight	378.32
Soluble In	Water
Chemical Structure	



Spectral Properties

Excitation Wavelength	571 nm
Emission Wavelength	584 nm

Applications

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 in sources that do not also contain high levels of phospholipase C.