

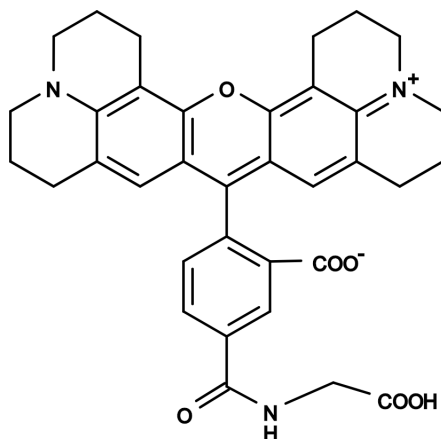
## Product Information Sheet

### Ordering Information

Product Number:	396
Product Name:	5-ROX glycine *Fluorescence reference standard for PCR reactions*
Unit Size:	25 mg
Storage Conditions:	Freeze (<-15 °C), Desiccated, Avoid Light
Expiration Date:	12 months upon receiving

### Chemical and Spectral Properties

Appearance:	Solid
Molecular Weight:	591.65
Chemical Structure:	



Soluble In:	Water
Excitation Wavelength:	575
Emission Wavelength:	602

### Application Notes

5-ROX glycine is used as a reference dye for performing PCR detections. Compared to the standard ROX dyes 5-ROX glycine has improved stability. The 5-ROX glycine has the similar spectral properties to those of 6-ROX.