

## 6-ROX glycine \*25 uM fluorescence reference solution for PCR reactions\*

Catalog Number: 395

Unit Size: 5 ml

### Product Details

---

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

### Chemical Properties

---

|                    |  |
|--------------------|--|
| Appearance         | Liquid dark red  |
| Molecular Weight   | 591.65   |
| Soluble In         | Water  |
| Chemical Structure |  |

### Spectral Properties

---

|                       |        |
|-----------------------|--------|
| Excitation Wavelength | 578 nm |
| Emission Wavelength   | 604 nm |

### Applications

---

6-ROX is predominately used as a reference dye for performing PCR detections. However, 6-ROX is very unstable compared to other rhodamine dyes. 6-ROX glycine has improved stability. The 6-ROX glycine has the similar spectral properties to those of 6-ROX. This 6-ROX dye is 25 uM solution in 20 mM Tris (pH 8.4), 0.1 mM EDTA and 0.01% Tween® 20.