

## Product Information Sheet

### Ordering Information

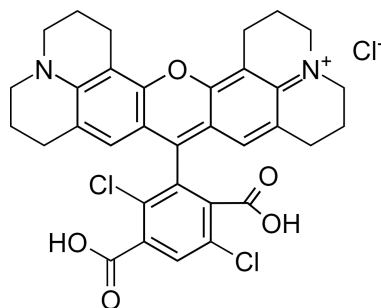
Product Number:	311
Product Name:	6-dROX [6-Carboxy-4,7-dichloror-X-hodamine]
Unit Size:	25 mg
Storage Conditions:	Freeze (<-15 °C), Desiccated, Avoid Light
Expiration Date:	12 months upon receiving

### Chemical and Spectral Properties

Appearance: Solid

Molecular Weight: 639.95

Chemical Structure:



Soluble In: DMSO

Excitation Wavelength: 602

Emission Wavelength: 621

### Application Notes

6-dROX (6-Carboxy-4,7-dichloror-X-hodamine) is one of the four critical 4,7-dichloro-rhodamine dyes used as the BigDye terminators for DNA sequencing in combination with dR110, dR6G and dROX. ABI PRISM® BigDye™ Terminator Cycle Sequencing is the classic method for DNA sequencing.