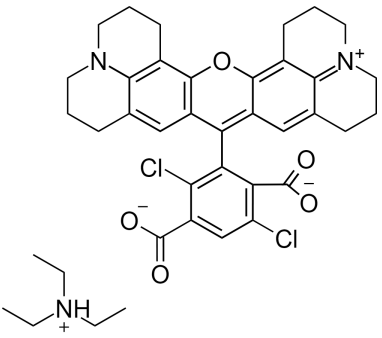
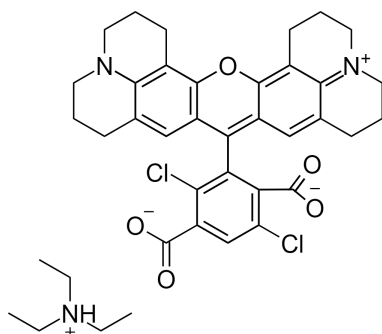


6-dROX [6-Carboxy-4,7-dichloror-X-hodamine]Catalog number: 311
Unit size: 25 mg**Product Details**

Storage Conditions	Freeze ($-15\text{ }^{\circ}\text{C}$), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	704.69
Soluble In	DMSO
Chemical Structure	

**Applications**

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.