

Product Information Sheet

Ordering Information

Product Number: 300

Product Name: 5-dR6G [5-Carboxy-4,7-dichlororhodamine 6G]

Unit Size: 25 mg

Storage Conditions: Freeze (<-15 °C), Desiccated, Avoid Light

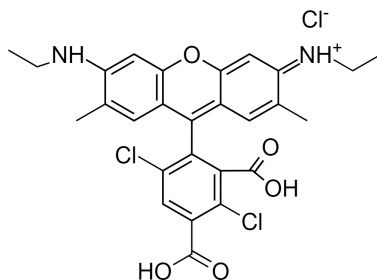
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Appearance: Solid

Molecular Weight: 563.86

Chemical Structure:



Soluble In: DMSO

Excitation Wavelength: 548

Emission Wavelength: 567

Application Notes

5-dR6G (5-Carboxy-4,7-dichlororhodamine 6G) is one of the four critical 4,7-dichloro-rhodamine dyes used as the BigDye terminators for DNA sequencing in combination with dR110, dTMR and dROX. ABI PRISM® BigDye™ Terminator Cycle Sequencing is the classic method for DNA sequencing.