

Catalog #:AS-81106

AnaSpec, Inc.

Analytical Data Sheet

Product Name:	6-CR 6G,SE (6-carboxyrhodamine 6G, succinimidyl ester)
Lot No.:	CM 9040
Weight:	5mg
Formula:	$C_{31}H_{29}N_3O_7$
Molecular Weight:	555.59
Appearance:	Dark red solid
Purity:	
TLC:	>90%
HPLC:	>90% purity

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ **Date:** _____

Quality Control