

Catalog #: 81105

AnaSpec, Inc.

Analytical Data Sheet

Product Name: 5-CR 6G, SE(5-Carboxyrhodamine 6G, succinimdy ester

Lot No.: CM5120

Weight: 5 mg

Formula: $C_{31}H_{29}N_3O_7$

Molecular Weight: 555.59

Appearance: Dark red solid

Purity:

HPLC: >90 %

TLC: >90%

Remarks: *Not for Human Use. Research Purposes Only.*

Release By: _____ **Date:** _____

Quality Control