

Catalog #:AS-81130

AnaSpec, Inc.

Analytical Data Sheet

Product Name: Sulforhodamine 101 sulfonyl chloride

Lot No.: CM2081

Weight: 10 mg

Appearance: Dark purple solid

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

Released By: _____ **Date:** _____
Quality Control