

Catalog #: 81833

AnaSpec, Inc.

Analytical Data Sheet

Product Name: QXL™ 520 C2 maleimide, disodium salt

Lot No.:

Weight: 10mg

Formula: N/A

Molecular Weight: 675.60

Appearance: Red solid

Purity:

HPLC: 90%

TLC: >90%

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

Release By: _____ **Date:** _____

Quality Control