

Catalog #: 81828

AnaSpec, Inc.

Analytical Data Sheet

Product Name: QXL™490 C₂ amine

Lot No.:

Weight: 10 mg

Formula:

Molecular Weight: 533.53

Appearance: Red solid

Purity:

TLC: >95%

HPLC: 95%

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ **Date:** _____

Quality Control