

Catalog #: 81823

AnaSpec, Inc.

Analytical Data Sheet

Product Name: DABCYL Plus™ C₂ amine

Lot No.:

Weight: 10 mg

Appearance: Red solid

Ninhydrin Test:

Substitution Level:

Particle Size:

Pore Size:

Structure:

Additional Instructions: For best results, pre-swell resin with DCM for 10 minutes.

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ **Date:** _____
Quality Control