

Catalog #: 81121

AnaSpec, Inc.

Analytical Data Sheet

Product Name: 5-TAMRA(5-carboxytetramethyl rhodamine)

Lot No.: CM 3012

Weight: 10 mg

Appearance: Dark 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