

Catalog #: 29284

AnaSpec, Inc.

Analytical Data Sheet

Product Name: (S)-N-1-Boc-2-isopropylpiperazine (Stock#: C-1132)

Lot No.: CE10645

Weight: 5 g

Appearance: Liquid

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