

Catalog #: 26283-B01

AnaSpec, Inc.

Analytical Data Sheet

Product Name: Boc- α -methyl-L-aspartic acid

Lot No.: n/a

Weight: 100 mg

Formula: C₁₀H₁₇NO₆

Molecular Weight: 247.3

Appearance: n/a

Storage: Keep in a cool & dry place

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

Release By: _____ **Date:** _____

Quality Control