

Catalog #: 26432-B5

AnaSpec, Inc.

Analytical Data Sheet

Product Name: Boc- α -methyl-D-phenylalanine

Lot No.: AA32265

Weight: 5 g

Formula: C₁₅H₂₁NO₄

Molecular Weight: 279.3

Appearance: White crystalline powder

HPLC: $\geq 98\%$

Solubility: 1mmol in 1ml DMF

Storage: Keep in a cool & dry place

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

Release By: _____ **Date:** _____

Quality Control