

Catalog #:AS-29953-F1

AnaSpec, Inc.

Analytical Data Sheet

Product Name: Fmoc-N-Me-Ser(tBu)-OH

Lot No.: BG6106

Weight: 1 g

Formula: C₂₃H₂₇NO₅

Molecular Weight: 397.5

Appearance: Off-white solid

HPLC: ≥95%

Storage: Keep in a cool & dry place

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

Release By: _____ **Date:** _____

Quality Control