

Catalog #: 23491

AnaSpec, Inc.

Analytical Data Sheet

Product Name: Boc-Lys(DABCYL)-OH

Lot No.: AA34403

Weight: 5 g

Formula: $C_{26}H_{35}N_5O_5$

Molecular Weight: 497.6

Appearance: Yellow Solid

HPLC: 96%

Solubility: 1 mmol in 1 ml DMF

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

Release By: _____ **Date:** _____

Quality Control