

Catalog #: 23496

AnaSpec, Inc.

Analytical Data Sheet

Product Name:	Fmoc-Lys(DABCYL)-OH
Lot No.:	AA25399
Weight:	1 g
Formula:	$C_{36}H_{37}N_5O_5$
Molecular Weight:	619.7
Appearance:	Orange powder
HPLC:	≥95%
Solubility:	1mmol in 1ml DMF
Storage:	Store in a cool , dry & dark place

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ **Date:** _____

Quality Control