

Catalog #: 26274-F5

AnaSpec, Inc.

Analytical Data Sheet

Product Name:	Fmoc-L-2-aminoadipic acid
Lot No.:	AA22031
Weight:	5 g
Formula:	$C_{21}H_{21}NO_6$
Molecular Weight:	383.5
Appearance:	white powder
Purity:	
HPLC:	≥95% purity
TLC:	one spot (DCM:MeOH:AcOH = 9:1:0.1)
Solubility:	1mmol in 1ml DMF
Storage:	keep in a cool and dry place

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ **Date:** _____

Quality Control