

Catalog #: 23497

AnaSpec, Inc.

Analytical Data Sheet

Product Name: Fmoc-Lys(DABCYL)-OH

Lot No.: AA12886

Weight: 5 g

Formula: $C_{36}H_{37}N_5O_5$

Molecular Weight: 619.7 (Found : 620.9 / M + H⁺)

Appearance: Red Powder

Purity:

HPLC: > 95 %

TLC: one spot (CHCl₃ : MeOH = 9 : 1)

Solubility: 1 mmol in 10 ml DMF

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ **Date:** _____

Quality Control