

Catalog #: 23496

# AnaSpec, Inc.

## Analytical Data Sheet

---

**Product Name:** Fmoc-Lys(DABCYL)-OH

**Lot No.:** AA12886

**Weight:** 1 g

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

**Molecular Weight:** 619.7 (Found : 620.9 / M + H<sup>+</sup>)

**Appearance:** Red Powder

**Purity:**

**HPLC:** > 95 %

**TLC:** one spot (CHCl<sub>3</sub> : MeOH = 9 : 1)

**Solubility:** 1 mmol in 10 ml DMF

---

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

**Release By:** \_\_\_\_\_ **Date:** \_\_\_\_\_

*Quality Control*