

Catalog #: 23494

# AnaSpec, Inc.

## Analytical Data Sheet

---

**Product Name:** Fmoc-Glu(EDANS)-OH

**Lot No.:** AA12638

**Weight:** 1 g

**Formula:**  $C_{32}H_{31}N_3O_8S$

**Molecular Weight:** 617.7

**Appearance:** off-white powder

**HPLC:** >95%

**TLC:** one spot  $CHCl_3$ :MeOH 9:1

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

---

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

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

*Quality Control*