

Catalog #: 23489

**AnaSpec, Inc.**

## Analytical Data Sheet

---

**Product Name:** Boc-L-Glu(EDANS)-OH

**Lot No.:** AA8548

**Weight:** 5 g

**Formula:**  $C_{22}H_{29}N_3O_8S$

**Molecular Weight:** 495.6

**Appearance:** Off-white solid

**HPLC:**  $\geq 90\%$

**Solubility:** dissolve 1 mg of sample in 1.0 mL DMF .

**Storage:** keep in a dry & cool place .

---

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

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

*Quality Control*