

Catalog #: 26371-F05

**AnaSpec, Inc.**

## Analytical Data Sheet

---

**Product Name:** Fmoc-3-exo-aminobicyclo(2.2.1)hept-5-ene-2-exo-carboxylic acid

**Lot No.:** AA33495

**Weight:** 0.5 g

**Formula:**  $C_{23}H_{21}NO_4$

**Molecular Weight:** 375.4

**Appearance:** White powder

**Purity:**  
**HPLC:**  $\geq 98\%$

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

**Storage:** Keep in a dark cool & dry place

---

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

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

*Quality Control*