

Catalog #:AS-26260-F1

AnaSpec, Inc.

Analytical Data Sheet

---

**Product Name:** Fmoc-tert-butyl-L-alanine

**Lot No.:** AA32175

**Weight:** 1 g

**Formula:** C<sub>22</sub>H<sub>25</sub>NO<sub>4</sub>

**Molecular Weight:** 367.4

**Appearance:** White powder

**Purity:**

**HPLC:** 95%

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

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

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

---

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

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

*Quality Control*