

Catalog #: 26046-B5

**AnaSpec, Inc.**

## Analytical Data Sheet

---

**Product Name:** Boc-3-amino-2-methylbenzoic acid

**Lot No.:** AA33637

**Weight:** 5 g

**Formula:**  $C_{13}H_{17}NO_4$

**Molecular Weight:** 251.3

**Appearance:** Light purple powder

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

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

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

---

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

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

*Quality Control*