

Catalog #: 29272

# AnaSpec, Inc.

## Analytical Data Sheet

---

**Product Name:** (+)-N-4-Boc-2-methylpiperazine

**Lot No.:** CE11165

**Weight:** 5 g

**Formula:**  $C_{10}H_{20}N_2O_2$

**Molecular Weight:** 200.3

**Appearance:** Liquid

**Purity:**

**HPLC:** 98 %

**\* Absorbs CO<sub>2</sub> from air!**

**Handle and Store:** under nitrogen!

---

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

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

*Quality Control*