

Catalog #: 26247-B1

**AnaSpec, Inc.**

## Analytical Data Sheet

---

**Product Name:** Boc-D-cyclohexylglycine-OH · H<sub>2</sub>O

**Lot No.:** AG0641

**Weight:** 1 g

**Formula:** C<sub>13</sub>H<sub>23</sub>NO<sub>4</sub> · H<sub>2</sub>O

**Molecular Weight:** 257.3 · 18.0

**Appearance:** White powder

**Optical Rotation:**  $[\alpha]_D^{20} = -12.6^\circ$  (C=1 in MeOH corrected for H<sub>2</sub>O)

**TLC:** ≥ 99 %

**Storage:** Store at 2-8°C

---

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

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

*Quality Control*