

Catalog #: AS-21304 (1 g);  
AS-21304-5 (5 g)

AnaSpec, Inc.

## Analytical Data Sheet

---

Product Name:	Boc-4-Chloro-Phe-OH
Lot No.:	BG0210
Formula:	$C_{14}H_{18}NO_4Cl$
Molecular Weight:	299.8
Appearance:	Off-white powder
Optical Rotation:	$[\alpha]_D^{20} = +14^{\circ}C$ (C = 1, MeOH)
HPLC:	$\geq 98\%$
Storage:	Keep in a dry cool place

---

Remarks: Not for Human Use. Research Purposes Only.

Released By: \_\_\_\_\_ Date: \_\_\_\_\_  
Quality Control