

Catalog #: 26341-B1

AnaSpec, Inc.

Analytical Data Sheet

Product Name: Boc-(±)-4-amino 3-hydroxybutanoic acid

Lot No.: AA20168

Weight: 1 g

Formula: C₉H₁₈NO₅

Molecular Weight: 219.1

Appearance: white powder

HPLC: ≥ 98% purity

TLC: one spot (DCM:MeOH:AcOH 9:1:0.1)

Solubility: 1 mmol in 1ml DMF

Storage: Store in a cool dry place

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

Release By: _____ **Date:** _____

Quality Control