

Catalog #: 26197-B5

AnaSpec, Inc.

Analytical Data Sheet

Product Name: Boc-4-carboxymethylpiperazine

Lot No.: AA10901

Weight: 5 g

Formula: $C_{11}H_{20}N_2O_4$

Molecular Weight: 244.3

Appearance: off-white powder

Melting Point: 244 - 245°C

HPLC: $\geq 96\%$

Loss on Drying: n.d.

Free Amino Acids: $\leq 0.1\%$

Storage: keep dark, cool and dry

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

Release By: _____ **Date:** _____

Quality Control