

Catalog #: 26197-F1

AnaSpec, Inc.

Analytical Data Sheet

Product Name: Fmoc-4-carboxymethylpiperazine

Lot No.: AA25169

Weight: 1 g

Formula: $C_{21}H_{22}N_2O_4$

Molecular Weight: 366.4

Appearance: white powder

HPLC: $\geq 95\%$

Solubility: 1 mmol in 1 mL DMF.

Storage: Keep in a dry cool place.

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ **Date:** _____

Quality Control