

Catalog #: 81811

AnaSpec, Inc.

Analytical Data Sheet

Product Name: DABSYL-L-Methionine

Lot No.:

Weight: 25mg

Formula: $C_{19}H_{24}N_4O_4S_2$

Molecular Weight: 436.55

Appearance: Orange Solid

Purity:

HPLC: >95%

TLC: >95%

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ **Date:** _____

Quality Control