

Catalog #: 81813

AnaSpec, Inc.

Analytical Data Sheet

Product Name: DABSYL-L-Tryptophan

Lot No.:

Weight: 25mg

Formula: $C_{25}H_{25}N_5O_4S$

Molecular Weight: 491.56

Appearance: Orange Solid

Purity:

HPLC: >95%

TLC: >95%

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ **Date:** _____

Quality Control