

Catalog #: 23488

AnaSpec, Inc.

Analytical Data Sheet

Product Name: Boc-L-Glu(EDANS)-OH

Lot No.: AA8548

Weight: 1 g

Formula: $C_{22}H_{29}N_3O_8S$

Molecular Weight: 495.6

Appearance: Off-white solid

HPLC: ≥ 90 %

Solubility: dissolve 1 mg of sample in 1.0 mL DMF .

Storage: keep in a dry & cool place .

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ **Date:** _____

Quality Control