

Catalog #: 23486-01

AnaSpec, Inc.

Analytical Data Sheet

Product Name: Boc-Asp(EDANS)-OH

Lot No.: AA29279

Weight: 0.1 g

Formula: $C_{21}H_{27}N_3O_8S$

Molecular Weight: 481.6

Appearance: Light yellow powder

HPLC: $\geq 85\%$

Solubility: 1mmol in 1ml DMF

Storage: Keep in a dry cool place .

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

Release By: _____ **Date:** _____

Quality Control