

Catalog #: 23492-01

AnaSpec, Inc.

Analytical Data Sheet

Product Name: Fmoc-Asp(EDANS)-OH

Lot No.: AA 19760

Weight: 100 mg

Formula: $C_{31}H_{29}N_3O_8S$

Molecular Weight: 603.7

Appearance: Grey Powder

HPLC: $\geq 95\%$ purity

TLC: One spot (CHCl₃:MeOH=9:1)

Solubility: 1 mmol in 10ml DMF

Storage: Store in a cool & dry place

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

Release By: _____ **Date:** _____

Quality Control