

Catalog #: 23487

AnaSpec, Inc.

Analytical Data Sheet

Product Name:	Boc-Asp(EDANS)-OH
Lot No.:	AA29279
Weight:	5 g
Formula:	$C_{21}H_{27}N_3O_8S$
Molecular Weight:	481.6
Appearance:	Light yellow powder
HPLC:	≥ 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