

Catalog #: AS-26172-B01

AnaSpec, Inc.

Analytical Data Sheet

Product Name:	Boc-6-fluoro-D,L-tryptophan
Lot No.:	AA25163
Weight:	0.1 g
Formula:	$C_{16}H_{19}FN_2O_4$
Molecular Weight:	322.3
Appearance:	White powder
Purity:	$\geq 98\%$ (HPLC)
TLC:	One spot (CHCl ₃ :MeOH:HOAc = 9:1:0.1)
Solubility:	1 mmol in 1 ml DMF
Storage:	Keep in a dark, cool and dry place

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ Date: _____
Quality Control