

Catalog #: 27131

AnaSpec, Inc.

Analytical Data Sheet

Product Name:	Fmoc-D-Arg(Pbf)-OH N- α -Fmoc-N ⁶ -2,2,4,6,7-pentamethyldihydrobenzofuran-5-sulfonyl-D-arginine
Lot No.:	BO13.1
Weight:	648.8 mg (1mmol)
Formula:	C ₃₄ H ₄₀ N ₄ O ₇ S
Molecular Weight:	648.8
Appearance:	white to off-white powder
Optical Rotation:	$[\alpha]_D^{25} = + 4.7^\circ$ (c=1, DMF)
HPLC:	$\geq 99\%$ purity
Diisopropylether content:	5.54 %
Water content:	0.97 %
Solubility:	Clear colorless solution in 5% m/v in DMF
Storage:	Store in a dry cool place

Remarks: Not for Human Use. Research Purposes Only.

Release By: _____ **Date:** _____

Quality Control