

(2R,3R,4R,5R)-2-((bis(4-methoxyphenyl)(phenyl)methoxy)methyl)-4-((tert-butyldimethylsilyl)oxy)-5-(5-methoxy-2,4-dioxo-3,4-dihydropyrimidin-1(2H)-yl)tetrahydrofuran-3-yl (2-cyanoethyl) diisopropylphosphoramidite

Purity Specification:	≥95%		
Molecular formula:	C46H63N4O10PSi		
CAS number:	Formula weight:	891.09	Expiration date: 09JAN2030

Tests	Results
HPLC	Purity: 95.4 %
Mass Spec	M-H: 889.6
NMR	Conforms
31P-NMR	Conforms

Gibby Perry

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM