

1-[(4R,5S)-5-[(2-aminoethyl-1,1,2,2-d4)amino]-N2-(10,12-dimethyl-1-oxotetradecyl)-4-hydroxy-L-ornithine]-5-[(3R)-3-hydroxy-L-ornithine]-pneumocandin B0, diacetate

Expiration date: 25SEP2029

Tests	Results
Elemental analysis	C: 50.79% H: 7.94% N: 10.34%
HPLC	Purity: 98.2 %
NMR	Conforms

Subburay

WWW.CAYMANCHEM.COM