

Cayman
CHEMICAL

(2R,2'R)-2,2'-((((5R,6R,7S,9S,11R,18S,19S)-19-amino-11,18-dihydroxy-5,9-di(methyl-13C)icosane-6,7-diyl-1,2,3,4,5,6,7,8,9,10,11,12,13,14,15,16,17,18,19,20-13C20)bis(oxy))bis(2-oxoethane-2,1-diyl-1,2-13C2))disuccinic-13C4 acid
*Item No. 31274 * Batch No. 0818437*

Purity Specification:	≥95%			
Molecular formula:	[13C]34H59NO14			
CAS number:	Formula weight:	739.60	Expiration date:	19MAR2030

Tests	Results
HPLC	Purity: 98.8 %
Concentration	10.00 ug/mL

Reviewed and approved by: Justine Albertson

Subburay

WWW.CAYMANCHEM.COM