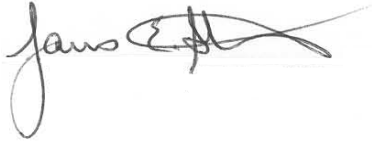


Certificate of Analysis

Catalogue number:	UM304	
Product name:	Biotin K63-TUBE1	
Lot number:	TR-41858.001	
Expiration Date:	One year after date of Purchase	
Package size:	50µg	
Form:	Liquid, 0.5mg/mL	
Buffer:	PBS	
Storage:	-80°C. Avoid repeated freeze/thaw cycles	
QC Manager:		
Protein:	Expected	Determined
Molecular weight	10,320.3 Da	10,321.3 Da
Method		MS
Specific Activity		
Method		
Purity (%)	>90	>98.8
Method		RP-HPLC
Vector:		
Form (circular/linear)		
Parent		
Size		
Sequence		N/A
Digest test		N/A
Insert test		N/A

Description:

Biotin K63-TUBE 1 is useful for detection and characterization of K63-linked polyubiquitylated proteins by Far Western blotting, affinity purification from lysates and biological fluids, and in situ detection.