


Certificate of Analysis

Catalogue number:	UM607
Product name:	K48 TUBE HF FLAG
Lot number:	UM-43138.001
Expiration Date:	1 year from date of receipt
Package size:	50µg, 250µg
Form:	Liquid
Buffer:	PBS, pH 7.2
Storage:	minus (-) 80 C. Avoid repeated freeze/thaw cycles
QC Manager:	

Protein:	Expected	Determined
Molecular weight	22994.21 Da	22996.16 Da
Method		MS
Specific Activity		n/a
Method		
Purity (%)	>90	94.65%
Method		RP-HPLC
Vector:		
Form (circular/linear)		
Parent		
Size		
Sequence		N/A
Digest test		N/A
Insert test		N/A