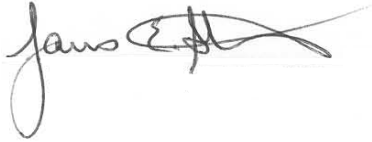


Certificate of Analysis

Catalogue number:	AB7030	
Product name:	Affinity purified, chicken anti-HuSUMO3 IgY	
Lot number:	AB-40911.001	
Expiration Date:	One year from date of purchase	
Package size:	50µg	
Form:	liquid, 0.5mg/mL	
Buffer:	Tris-buffered saline	
Storage:	-20°C or lower, avoid repeated freeze/thaw cycles	
QC Manager:		
Protein:	Expected	Determined
Molecular weight	not tested	
Method		MS
Reactivity	hSUMO3	10ng at 1:1000 dil'n
Method		Western Blot
Purity (%)	not tested	
Method		
Vector:		
Form (circular/linear)		
Parent		
Size		
Sequence		N/A
Digest test		N/A
Insert test		N/A

Description:

0