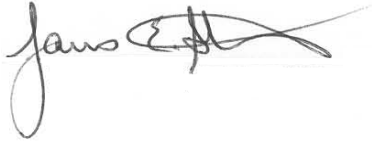


Certificate of Analysis

Catalogue number:	SI6312	
Product name:	DUB-resistant K63-linked diubiquitin (Ub2)	
Lot number:	Ub-41662.001	
Expiration Date:	One year from date of purchase	
Package size:	100 µg @ 1 mg/ml	
Form:	liquid	
Buffer:	20 mM Tris pH 7.5, 0.15 M NaCl, 1 mM EDTA	
Storage:	-80°C	
QC Manager:		
Protein:	Expected	Determined
Molecular weight	17,123.6 Da	17,124.3 Da
Method		MS
Specific Activity		N/A
Method		
Purity (%)	>90	>95
Method		RP-HPLC
Vector:		
Form (circular/linear)		
Parent		
Size		
Sequence		N/A
Digest test		N/A
Insert test		N/A

Description:

DUB-resistant, K63-linked Di-ubiquitin (Ub2) contains an amino acid substitution near the C-terminus of the distal ubiquitin that renders it resistant to deubiquitylases (DUBs). DUB-resistant ubiquitin chains are ideal reagents for structural or binding studies with deubiquitylases or other ubiquitin binding proteins. They can also be used as inhibitors or to determine the linkage-specificity of DUBs. The product contains a wild-type C-terminus on the proximal ubiquitin.