
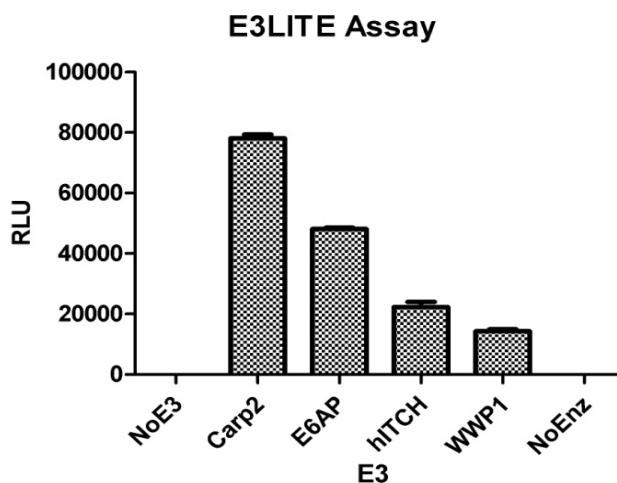
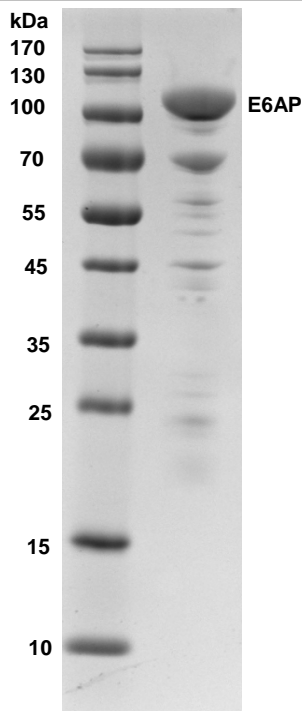


Certificate of Analysis

Catalogue number:	UB304T	
Product name:	His ₆ SUMO E6AP	
Lot number:	LG-41981.001	
Expiration Date:	One year from date of purchase	
Package size:	250µg	
Form:	Liquid, 17.5µM	
Buffer:	20mM Hepes pH7.5, 150mM NaCl, 10% Glycerol	
Storage:	-80°C	
QC Manager:		
Protein:	Expected	Determined
Molecular weight	112,838.5 Da	106.6kDa
Method		SDS-PAGE
Specific Activity		
Method		
Purity (%)		64.4%
Method		SDS-PAGE



all products are for research use only • not intended for human or animal diagnostic or therapeutic uses
 techsupport@lifesensors.com • www.lifesensors.com • sales@lifesensors.com
 Copyright © 2013 LifeSensors, Inc. All Rights Reserved