

K11-linked tri-ubiquitin (Ub3)

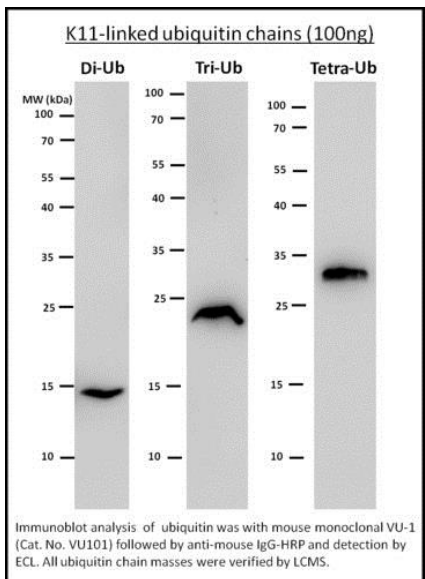
Cat. # SI1103

Background: Increasingly, researchers are focusing on the role poly-ubiquitin chains linked through K11. This post-translational modification has been linked to the ERAD cycle as a signal, similar to K48 linkage, for proteasomal degradation. More recently, K11 linkage appears to play an important role in cell cycle signaling, as it is associated with the anaphase promoting complex (APC) of ubiquitylation machinery.

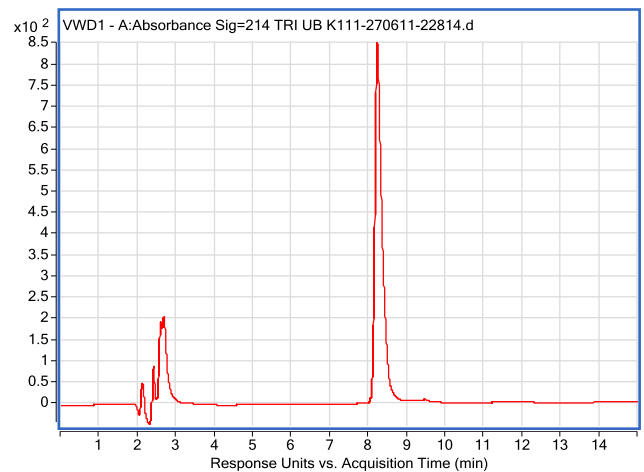
These tri-ubiquitin chains are generated from the enzymatic linkage (UBE2S) of wild-type ubiquitin through lysine 11 and purified to >95% homogeneity by ion exchange chromatography.

Product Information

Purity:	≥ 95% by Western blot analysis
Molecular Weight:	25,658.5 Da
Physical State:	Liquid; 20mM Tris-HCl, pH 7.5, 0.15M NaCl, 1mM EDTA
Quantity:	100 µg
Solubility:	>1mg/mL
Storage:	-80° C. Avoid repeated freeze/thaw cycles
Concentration:	Lot dependent, please see Certificate of Analysis or vial label for actual concentration



Western blot analysis



RP-HPLC Profile

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
 Copyright © 2009 LifeSensors, Inc. All Rights Reserved