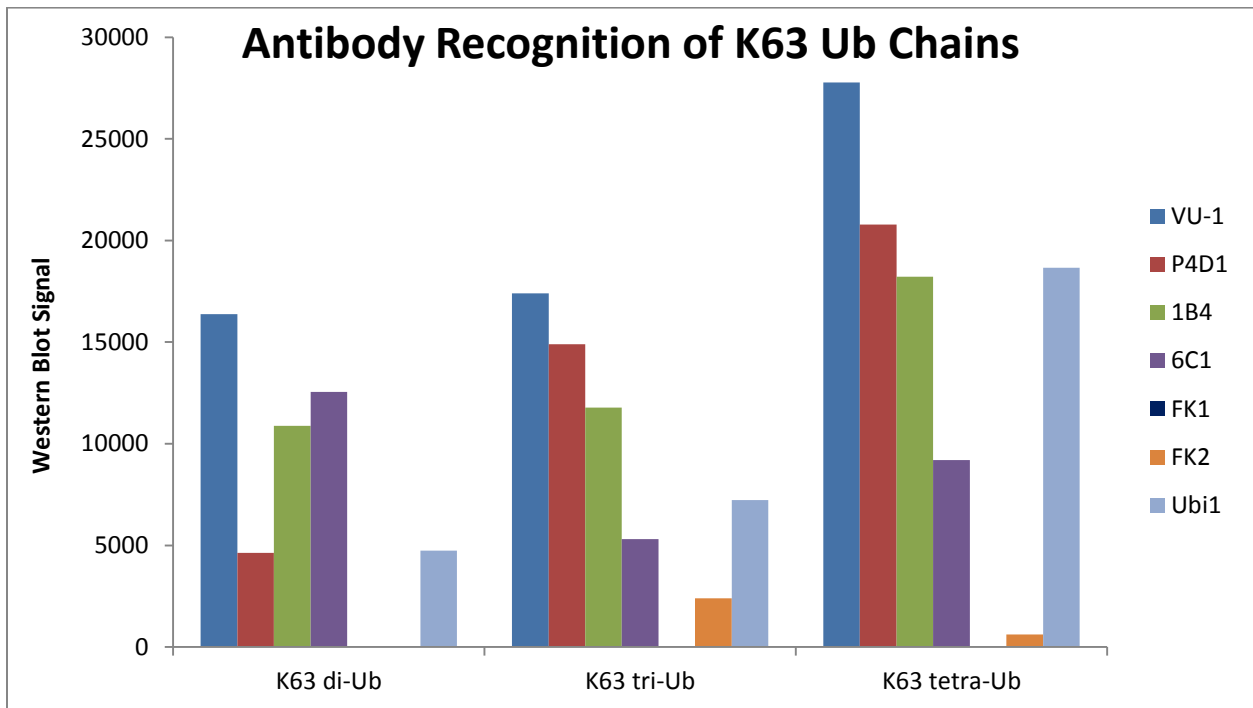
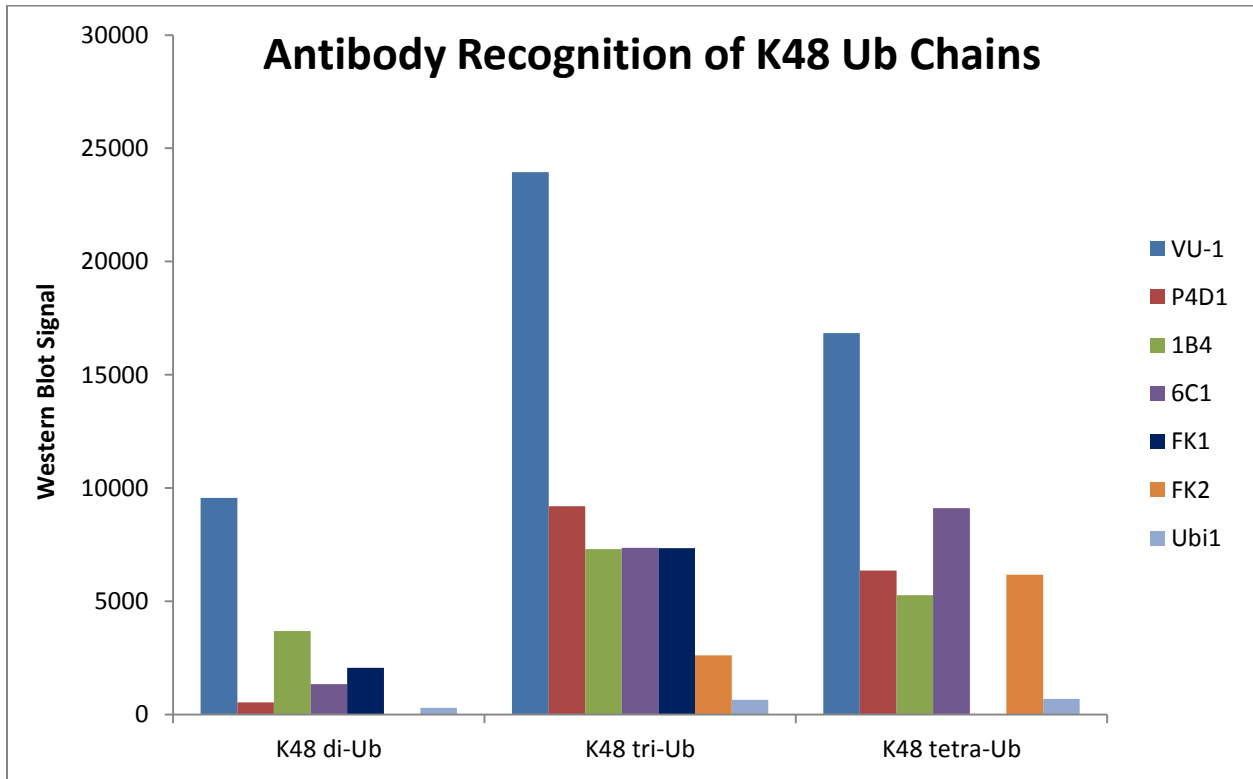
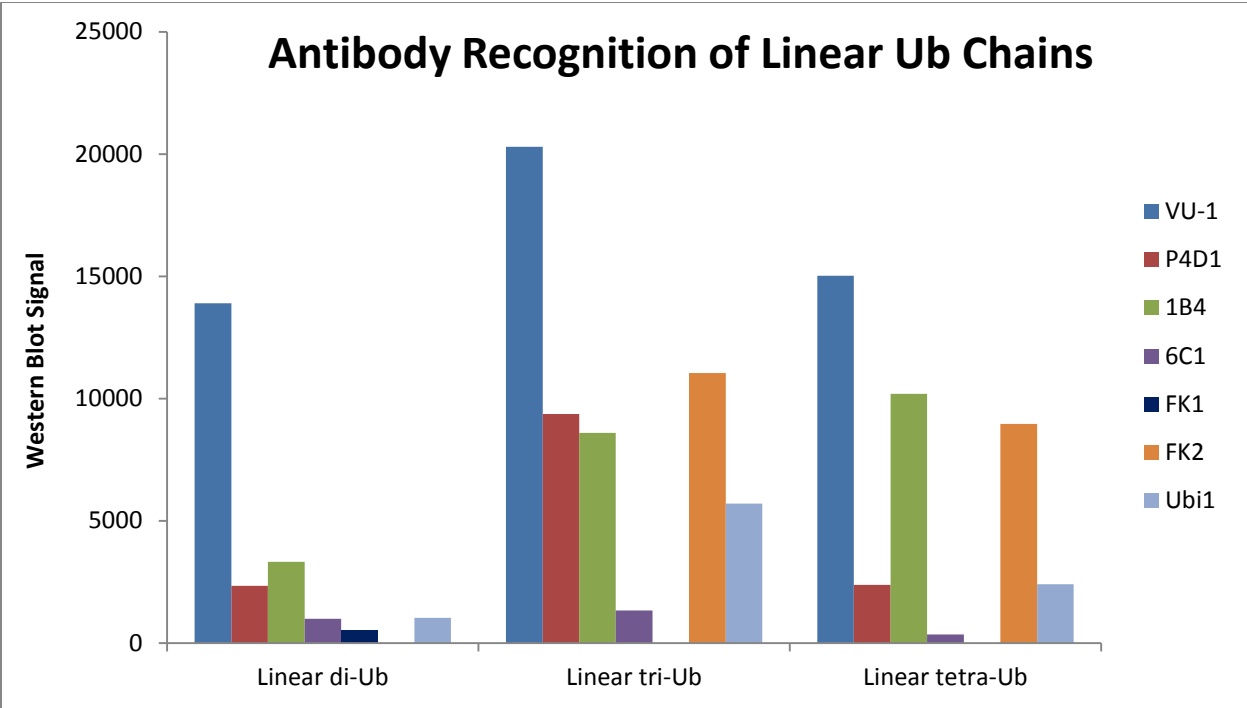
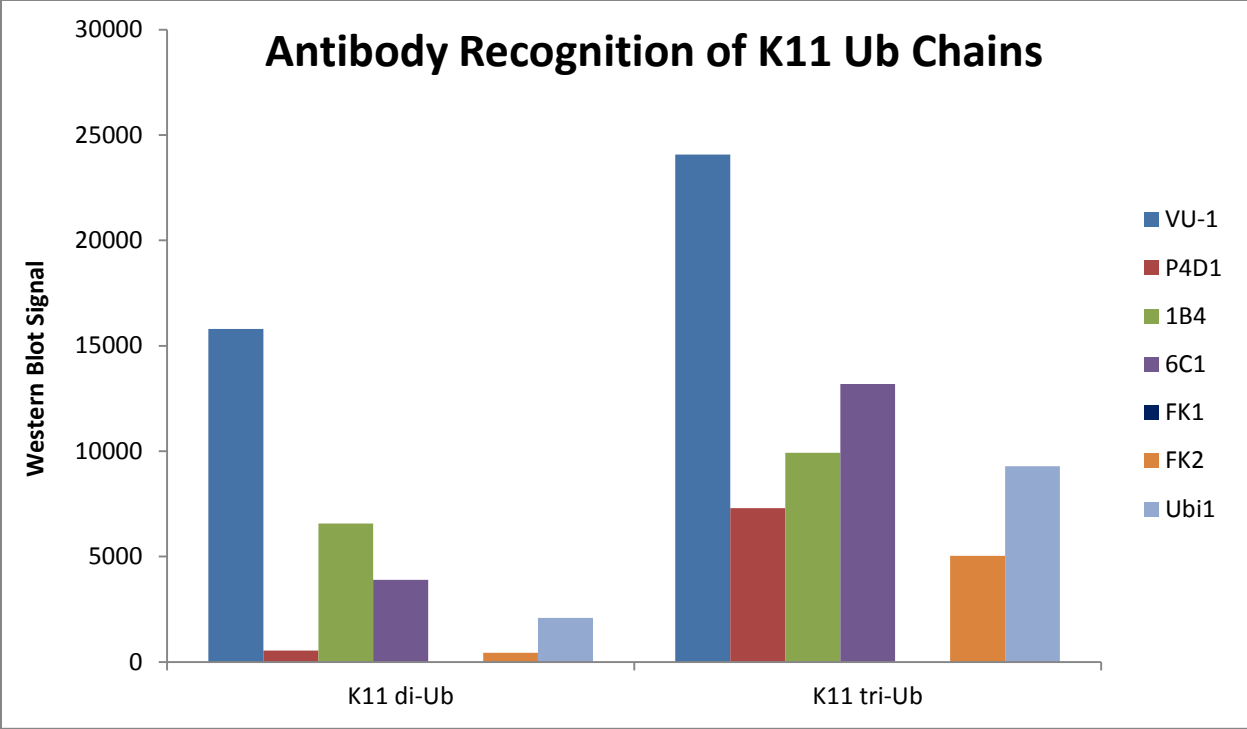


Table of Contents

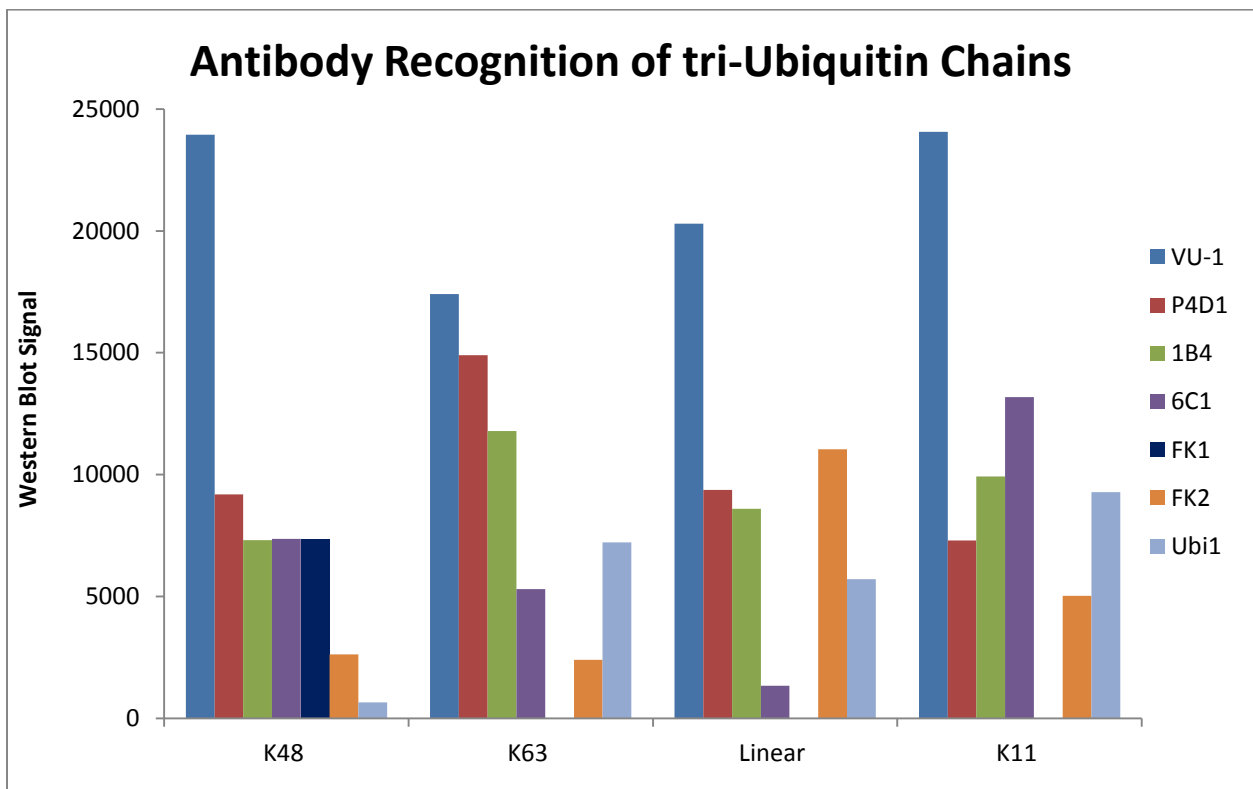
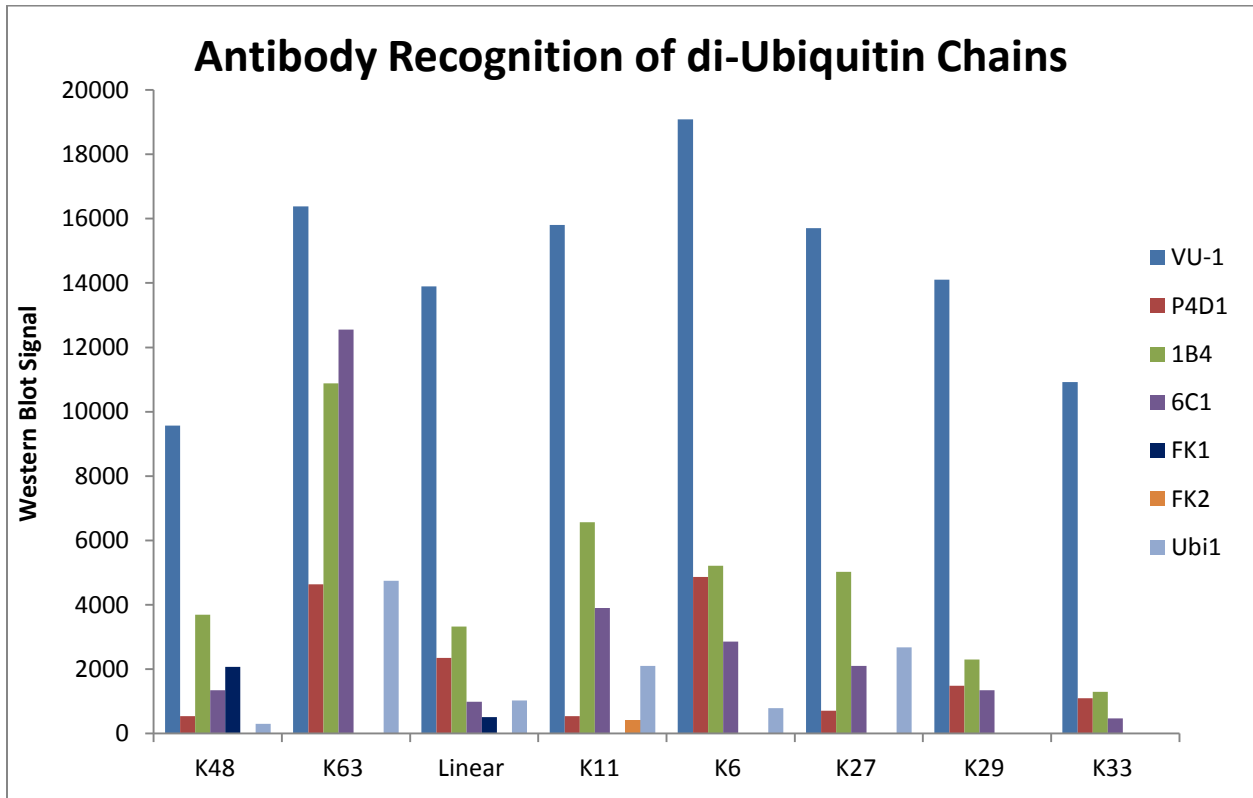
Antibody Comparison by Ub Chain Linkages.....	2
Antibody Comparison by Ub Chain Length.....	4
Western Blot.....	6

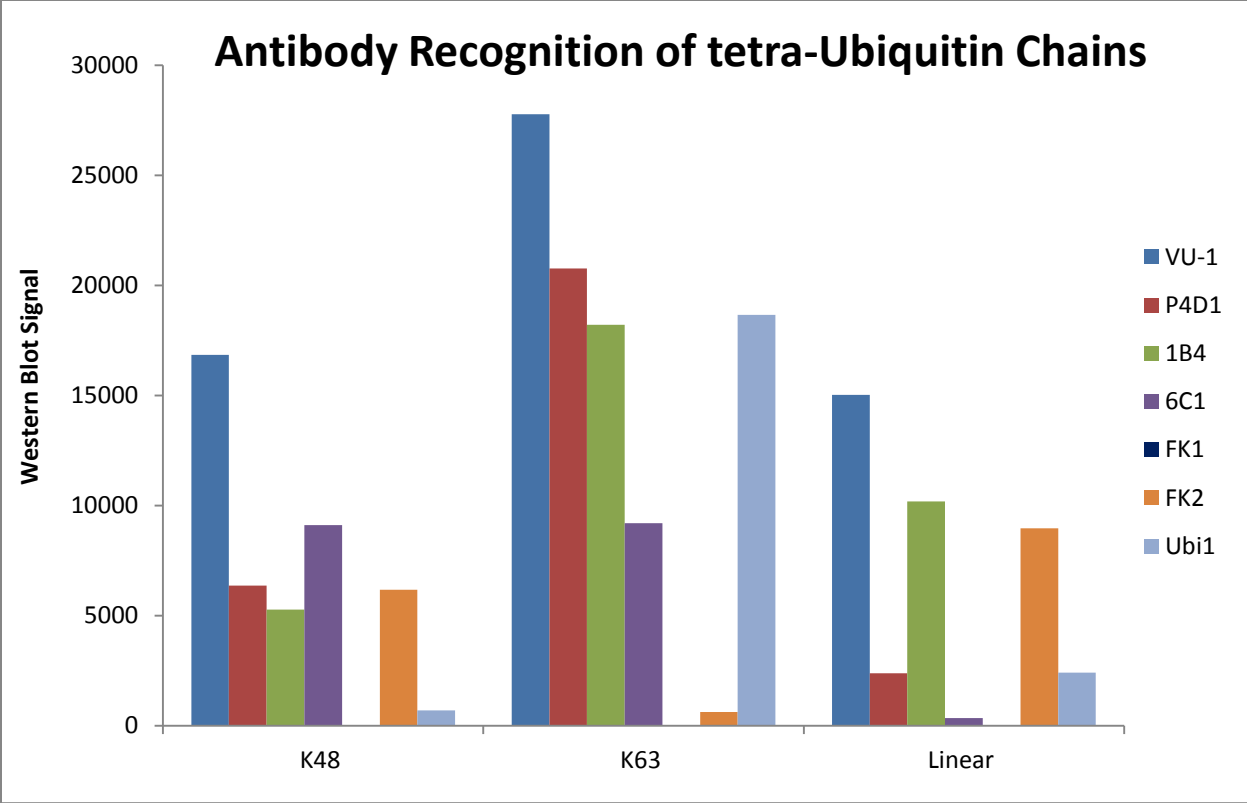
Antibody Comparison by Ub Chain Linkages





Antibody Comparison by Ub Chain Length





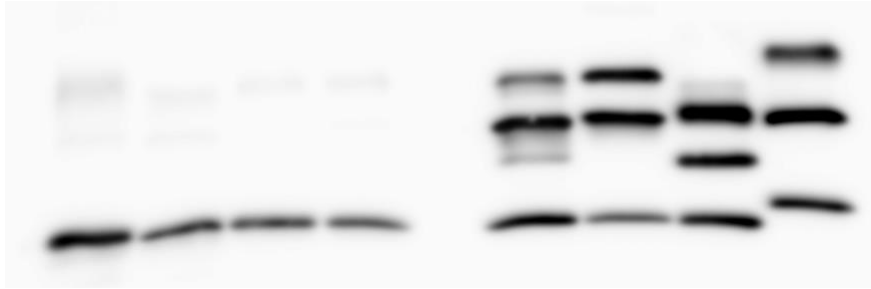
Western Blots

K6, K27, K29, K33: di-ubiquitin (20ng each)

K11: di- and tri-ubiquitin (20ng each)

K48, K63, linear: di-,tri-, and tetra-ubiquitin (20ng each)

K6 K27 K29 K33 K11 K48 K63 Linear



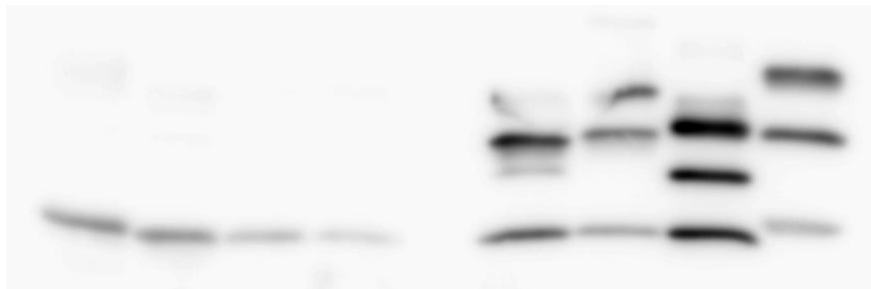
VU-1 antibody

K6 K27 K29 K33 K11 K48 K63 Linear



P4D1 antibody

K6 K27 K29 K33 K11 K48 K63 Linear



1B4 antibody

K6 K27 K29 K33 K11 K48 K63 Linear



6C1 antibody

K6 K27 K29 K33 K11 K48 K63 Linear



FK1 antibody

K6 K27 K29 K33 K11 K48 K63 Linear



FK2 antibody



Ubi1 antibody