

LifeSensors Products		Competitor Products			
		Millipore ®		Enzo Life Sciences ®	Cell Signaling®
Biotin k63-TUBE1		Anti-Ubiquitin, Lys63-Specific, clone Apu3 (05-1308)	Anti-Ubiquitin, Lys63-Specific, clone HWA4C4 (05-1313)	Polyubiquitin (K ⁶³ -linkage-specific), mAb (HWA4C4) (BML-PW0600)	K63-linkage Specific Polyubiquitin (D7A11) Rabbit mAb #5621
Quantity	200 µg	100 µl	100 µg	100 µl	100 µl
Price	\$250.00	\$369.00	\$329.00	\$312.00	\$269.00
Number of Blots	20	10 blots		10 blots	10 blots
Cost per Blot	\$12.50	\$36.90		\$31.20	\$26.90
Dilution	1 : 1000	1 : 1000	1 : 500	1 : 1000	Not Disclosed
Blot Denature/ Native	Native	Denature	Denature	Denature	Denature

¹ All prices and data are based on those published by companies as of 27th January 2012. Prices and information were retrieved from those published on websites.

² Hjerpe R, Aillet F, Lopitz-Otsoa F, Lang V, England P, Rodriguez MS., Efficient protection and isolation of ubiquitylated proteins using tandem ubiquitin-binding entities., *EMBO Rep.*, Pub Med: 2009 October 2

³ Assume 10ml per blot

Website Links

1) Millipore

- <http://www.millipore.com/catalogue/item/05-1308>
- <http://www.millipore.com/catalogue/item/05-1313>

2) Enzo Life Sciences

- <http://www.enzolifesciences.com/BML-PW0600/polyubiquitin-k63-linkage-specific-mab-hwa4c4/>

3) Cell Signaling

- <http://www.cellsignal.com/products/5621.html>

