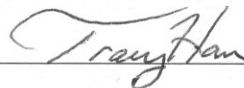


CERTIFICATE OF ANALYSIS

10X ELISA Wash Buffer (WB1)			
Catalog#	Size	Lot#	Expiry at room temperature
527	Bulk	14L45	07/2016
6251	25 mL	n/a	n/a
650	100 mL	n/a	n/a
651	500 mL	n/a	n/a
652	1 L	n/a	n/a
676	10 L	14L46	07/2016

Specifications	Criteria	Lot Data	Results
pH at 10X	6.8 – 7.0	6.80	Pass
pH at 1X	7.2 – 7.6	7.38	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	99,000 - 121,000 uMHO	107,633	Pass
Storage	Room temperature or refrigerated	Room temperature	Pass

Quality Control



Tracy Hanson

Date

07/29/2014

07/29/2014