

CERTIFICATE OF ANALYSIS

10X ELISA Wash Buffer (WB1)			
Catalog#	Size	Lot#	Expiry at room temperature
527	Bulk	14T48	10/2016
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	14T49	10/2016

Specifications	Criteria	Lot Data	Results
pH at 10X	6.8 – 7.0	6.81	Pass
pH at 1X	7.2 – 7.6	7.43	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	99,000 - 121,000 uMHO	105,633	Pass
Storage	Room temperature or refrigerated	Room Temperature	Pass

Quality Control



Tracy Hanson

Date

10/17/2014

10/17/2014