

# CERTIFICATE OF ANALYSIS

10X ELISA Wash Buffer (WB1)			
Catalog#	Size	Lot#	Expiry at room temperature
527	Bulk	15J53	01/2017
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	15J54	01/2017

Specifications	Criteria	Lot Data	Results
pH at 10X	6.8 – 7.0	6.81	Pass
pH at 1X	7.2 – 7.6	7.26	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity $\mu$ MHOs	99,000 - 121,000 $\mu$ MHO	109,500	Pass
Storage	Room temperature or refrigerated	Room Temperature	Pass

Quality Control



Tracy Hanson

Date

01/26/2015

01/26/2015