

CERTIFICATE OF ANALYSIS

10X ELISA Wash Buffer (WB1)			
Catalog#	Size	Lot#	Expiry at room temperature
527	Bulk	15G21	08/2017
6251	25 mL	NA	NA
650	100 mL	NA	NA
651	500 mL	NA	NA
652	1 L	NA	NA
676	10 L	15G22	08/2017

Specifications	Criteria	Lot Data	Results
pH at 10X	6.8 – 7.0	6.87	Pass
pH at 1X	7.2 – 7.6	7.45	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	99,000 - 121,000 uMHO	107,300	Pass
Storage	Room temperature or refrigerated	Room temperature	Pass

Quality Control

Tracy Hanson

Date

08/12/2015

Tracy Hanson

08/12/2015