

CERTIFICATE OF ANALYSIS

10X ELISA Wash Buffer (WB1)			
Catalog#	Size	Lot#	Expiry at room temperature
527	Bulk	18N49	06/2020
6251	25 mL	NA	NA
650	100 mL	NA	NA
651	500 mL	NA	NA
652	1 L	18N50	06/2020
676	10 L	NA	NA

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	115,367	Pass
Storage	Room temperature or refrigerated	Room temperature	Pass

Quality Control



Tracy Murphy

Date

06/29/2018

06/29/2018