

CERTIFICATE OF ANALYSIS

10X ELISA Wash Buffer (WB1)			
Catalog#	Size	Lot#	Expiry at room temperature
527	Bulk	18Y18	05/2020
6251	25 mL	NA	NA
650	100 mL	NA	NA
651	500 mL	NA	NA
652	1 L	18Y19	05/2020
676	10 L	18Y20	05/2020

Specifications	Criteria	Lot Data	Results
pH at 10X	6.8 – 7.0	6.88	Pass
pH at 1X	7.2 – 7.6	7.46	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMH0s	99,000 - 121,000 uMHO	114,800	Pass
Storage	Room temperature or refrigerated	Room temperature	Pass

Quality Control

Tracy Murphy

Tracy Murphy

Date

05/21/2018

05/21/2018