

CERTIFICATE OF ANALYSIS

10X ELISA Wash Buffer (WB1)			
Catalog#	Size	Lot#	Expiry at room temperature
527	Bulk	15B49	11/2017
6251	25 mL	NA	NA
650	100 mL	NA	NA
651	500 mL	NA	NA
652	1 L	NA	NA
676	10 L	15B50	11/2017

Specifications	Criteria	Lot Data	Results
pH at 10X	6.8 – 7.0	6.87	Approved
pH at 1X	7.2 – 7.6	7.47	Approved
Appearance	Clear liquid	Conforms	Approved
Conductivity uMHOs	99,000 - 121,000 uMHO	108,133	Approved
Storage	Room temperature or refrigerated	Refrigerated	Approved

Quality Control

Tracy Murphy

Tracy Murphy

Date

11/30/2015

11/30/2015