

CERTIFICATE OF ANALYSIS

Antigen Coating Buffer, 5X			
Catalog#	Size	Lot#	Expiry at Room Temperature
5136	Bulk	19T16	10/2021
6247	100 mL	19T17	10/2021
6248	500 mL	19T18	10/2021
6249	1 L	19T19	10/2021
6250	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH at 5X	7.8-8.2	8.13	Pass
pH at 1X	8.1-8.6	8.17	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity μ MHOs at 5X	45,000-55,000	51,100	Pass
Storage	Refrigerated or room temperature	Room temperature	Pass
Documentation	Approved	Approved	Pass

This is an assurance that the specified lot conforms to all requirements of the product specifications.
 This information is believed to be accurate and is offered in good faith.
 Product is for research use only. Not for use in diagnostic procedures.

Quality Assurance: Tracy Murphy
 Tracy Murphy

Date: 10/8/19
 10/08/2019