

CERTIFICATE OF ANALYSIS

Antibody Coating Buffer, 5X			
Catalog#	Size	Lot#	Expiry at Room Temperature
524	Bulk	19G27	08/2021
644	100 mL	NA	NA
645	500 mL	19G28	08/2021
646	1 L	19G29	08/2021
658	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH at 5X	7.8-8.2	8.17	Pass
pH at 1X	8.1-8.6	8.16	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity μ MHOs at 5X	45,000-55,000	50,400	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: 8/14/2019
 8/14/2019