

# CERTIFICATE OF ANALYSIS

<b>Neptune Blocking Buffer (BB2)</b>			
Catalog#	Size	Lot#	Expiry
53	Bulk	19G11	08/2021
62	100 mL	NA	NA
63	500 mL	19G12	08/2021
64	1 L	19G13	08/2021
660	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.41	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity uMH0s	15,000-19,000	16,350	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy  
 Tracy Murphy

Date 8/5/19  
 8/5/2019