

# CERTIFICATE OF ANALYSIS

## Neptune Blocking Buffer (BB2)

Catalog#	Size	Lot#	Expiry
53	Bulk	18J41	01/2020
62	100 mL	18J42	01/2020
63	500 mL	18J43	01/2020
64	1 L	18J44	01/2020
660	10 L	N/A	N/A

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.28	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity uMH0s	15,000-19,000	18,440	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

*Tracy Murphy*  
 Tracy Murphy

Date

*01/25/2018*  
 01/25/2018