

# CERTIFICATE OF ANALYSIS

## Neptune Blocking Buffer (BB2)

Catalog#	Size	Lot#	Expiry
53	Bulk	20J32	01/2022
62	100 mL	20A7	01/2022
63	500 mL	NA	NA
64	1 L	20J33	01/2022
660	10 L	N/A	N/A

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.49	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity uMHOs	15,000-19,000	16,720	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Sally Hed Dahlquist  
 Sally Hed Dahlquist

Date 4/15/2020  
 4/15/2020