

CERTIFICATE OF ANALYSIS

Neptune Blocking Buffer (BB2)

Catalog#	Size	Lot#	Expiry
53	Bulk	18D27	12/2020
62	100 mL	19Y39	12/2020
63	500 mL	NA	NA
64	1 L	18D28	12/2020
660	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.39	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity uMH0s	15,000-19,000	16,230	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

Tracy Murphy
 Tracy Murphy

Date

5/15/2019
 5/15/2019