

CERTIFICATE OF ANALYSIS

Neptune Blocking Buffer (BB2)

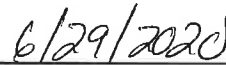
Catalog#	Size	Lot#	Expiry
53	Bulk	20N53	06/2022
62	100 mL	NA	NA
63	500 mL	NA	NA
64	1 L	20N54	06/2022
660	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.47	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity uMH0s	15,000-19,000	18,880	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


Sally Hed Dahlquist

Date


6/29/2020