

CERTIFICATE OF ANALYSIS

Neptune Blocking Buffer (BB2)


Catalog#	Size	Lot#	Expiry
53	Bulk	20Y27	05/2022
62	100 mL	20Y28	05/2022
63	500 mL	20Y29	05/2022
64	1 L	20Y30 20N51	05/2022 05/2022
660	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.44	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity uMH0s	15,000-19,000	16,290	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


 Sally Hed Dahlquist

Date


 6/26/2020