

CERTIFICATE OF ANALYSIS

Neptune Blocking Buffer (BB2)

Catalog#	Size	Lot#	Expiry
53	Bulk	18T6	10/2020
62	100 mL	NA	NA
63	500 mL	18T7	10/2020
64	1 L	NA	NA
660	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.37	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity uMHOs	15,000-19,000	16,050	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


 Kristi Strandberg

Date

10/8/18

10/8/2018