


# CERTIFICATE OF ANALYSIS

## General Blocking Buffer (BB1)

Catalog#	Size	Lot#	Expiry
517	Bulk	20J58	01/2022
632	100 mL	20J59	01/2022
633	500 mL	20J60	01/2022
640	1 L	20J61	01/2022
659	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.37	Pass
Appearance	Clear, slightly cloudy liquid	Conforms	Pass
Conductivity uMH0s	15,000-19,000	16,760	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

  
Tracy Murphy

Date

1/28/2020  
01/28/2020