

CERTIFICATE OF ANALYSIS

General Blocking Buffer (BB1)

Catalog#	Size	Lot#	Expiry
517	Bulk	20F24	02/2022
632	100 mL	NA	NA
633	500 mL	NA	NA
640	1 L	NA	NA
659	10 L	20F25	02/2022

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.35	Pass
Appearance	Clear, slightly cloudy liquid	Conforms	Pass
Conductivity uMHOs	15,000-19,000	16,730	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

Tracy Murphy
Tracy Murphy

Date

2/18/2020
2/18/2020