

CERTIFICATE OF ANALYSIS

General Blocking Buffer (BB1)


Catalog#	Size	Lot#	Expiry
517	Bulk	18S3	09/2020
632	100 mL	18S4	09/2020
633	500 mL	18S5	09/2020
640	1 L	18S6	09/2020
659	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.41	Pass
Appearance	Clear, slightly cloudy liquid	Conforms	Pass
Conductivity uMH0s	15,000-19,000	16,500	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


Kristi Strandberg

Date


09/04/2018