

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 20L14 20S10	01/2022 01/2022 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 uMHos	15,000-19,000	16,760	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Sally Hed Dahlquist
 Sally Hed Dahlquist

Date 9/4/2020
 9/04/2020