

CERTIFICATE OF ANALYSIS

General Blocking Buffer (BB1)			
Catalog#	Size	Lot#	Expiry
517	Bulk	20J44	01/2022
632	100 mL	NA	NA
633	500 mL	NA	NA
640	1 L	NA	NA
659	10 L	20J45	01/2022

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.38	Pass
Appearance	Clear, slightly cloudy liquid	Conforms	Pass
Conductivity uMH0s	15,000-19,000	17,020	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy
 Tracy Murphy

Date 1/22/2020
 1/22/2020