

CERTIFICATE OF ANALYSIS

General Blocking Buffer (BB1)

Catalog#	Size	Lot#	Expiry
517	Bulk	18D32	12/2020
632	100 mL	NA	NA
633	500 mL	NA	NA
640	1 L	NA	NA
659	10 L	18D33	12/2020

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.38	Pass
Appearance	Clear, slightly cloudy liquid	Conforms	Pass
Conductivity uMHOs	15,000-19,000	16,710	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy
 Tracy Murphy

Date 12/20/2018
 12/20/2018