

# CERTIFICATE OF ANALYSIS

## General Blocking Buffer (BB1)

Catalog#	Size	Lot#	Expiry
517	Bulk	16G69	08/2018
632	100 mL	NA	NA
633	500 mL	NA	NA
640	1 L	NA	NA
659	10 L	16G70	08/2018

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.41	Pass
Appearance	Clear, slightly cloudy liquid	Conforms	Pass
Conductivity uMHOs	15,000-19,000	17,180	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

*Tracy Murphy*  
Tracy Murphy

Date

*08/31/2016*  
08/31/2016