

# CERTIFICATE OF ANALYSIS

## General Blocking Buffer (BB1)

Catalog#	Size	Lot#	Expiry
517	Bulk	16L28	07/2018
632	100 mL	NA	NA
633	500 mL	16L31	07/2018
640	1 L	16L30	07/2018
659	10 L	16L29	07/2018

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.40	Pass
Appearance	Clear, slightly cloudy liquid	Conforms	Pass
Conductivity uMHOs	15,000-19,000	17,240	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

  
 Tracy Murphy

Date

07/20/2016  
 07/20/2016