



9401 James Avenue S., Suite #155, Bloomington, MN 55431 USA  
952-888-8788 fax 952-888-8988

Document: F11-517-4-B  
Effective: 03/19/13  
Supersedes: F11-517-4-A  
Page 1 of 1

# CERTIFICATE OF ANALYSIS

## General Blocking Buffer (BB1)

Catalog#	Size	Lot#	Expiry
517	Bulk	19Y32	05/2021
632	100 mL	19Y33 20J40	05/2021 05/2021
633	500 mL	19Y34	05/2021
640	1 L	19Y35	05/2021
659	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.36	Pass
Appearance	Clear, slightly cloudy liquid	Conforms	Pass
Conductivity uMHOs	15,000-19,000	16,690	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

  
Tracy Murphy

Date

1/16/2020  
1/16/2020