

CERTIFICATE OF ANALYSIS

General Blocking Buffer (BB1)

Catalog#	Size	Lot#	Expiry
517	Bulk	16R12	03/2018
632	100 mL	16R13	03/2018
633	500 mL	16R14	03/2018
640	1 L	16R15	03/2018
659	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.42	Pass
Appearance	Clear, slightly cloudy liquid	Conforms	Pass
Conductivity uMHOs	15,000-19,000	16,730	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


 Tracy Murphy

Date

03/09/2016
 03/09/2016