

CERTIFICATE OF ANALYSIS

General Blocking Buffer (BB1)			
Catalog#	Size	Lot#	Expiry
517	Bulk	16F41	02/2018
632	100 mL	NA	NA
633	500 mL	NA	NA
640	1 L	NA	NA
659	10 L	16F42	02/2018

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.40	Pass
Appearance	Clear, slightly cloudy liquid	Conforms	Pass
Conductivity uMH0s	15,000-19,000	15,680	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy
 Tracy Murphy

Date 02/18/2016
 02/18/2016