

# CERTIFICATE OF ANALYSIS

<b>SynBlock (BB3)</b>			
Catalog#	Size	Lot#	Expiry
512	Bulk	17R16	03/2019
641	100 mL	NA	NA
642	500 mL	NA	NA
643	1 L	NA	NA
661	10 L	17R17	03/2019

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.38	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	14,430	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control   
 Kristi Strandberg

Date 03/10/17  
 03/10/2017