

# CERTIFICATE OF ANALYSIS

<b>SynBlock (BB3)</b>			
Catalog#	Size	Lot#	Expiry
512	Bulk	18T8	10/2020
641	100 mL	NA	NA
642	500 mL	NA	NA
643	1 L	18T9	10/2020
661	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.30	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	14,220	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control   
 Kristi Strandberg

Date 10/19/18  
 10/9/2018