

# CERTIFICATE OF ANALYSIS

<b>SynBlock (BB3)</b>			
Catalog#	Size	Lot#	Expiry
512	Bulk	19R46	03/2021
641	100 mL	NA	NA
642	500 mL	NA	NA
643	1 L	NA	NA
661	10 L	19R47	03/2021

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.40	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	14,140	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy  
 Tracy Murphy

Date 3/26/19  
 3/26/2019