

# CERTIFICATE OF ANALYSIS

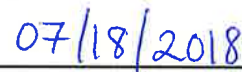
<b>SynBlock (BB3)</b>			
<b>Catalog#</b>	<b>Size</b>	<b>Lot#</b>	<b>Expiry</b>
512	Bulk	18Y14	07/2020
641	100 mL	NA	NA
642	500 mL	NA	NA
643	1 L	NA	NA
661	10 L	18Y15	07/2020

<b>Specifications</b>	<b>Criteria</b>	<b>Lot Data</b>	<b>Results</b>
pH	7.2-7.6	7.45	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	14,410	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

  
 Tracy Murphy

Date

  
 07/18/2018