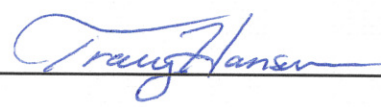


# CERTIFICATE OF ANALYSIS

<b>SynBlock (BB3)</b>			
<b>Catalog#</b>	<b>Size</b>	<b>Lot#</b>	<b>Expiry</b>
512	Bulk	14T7	10/2016
641	100 mL	N/A	N/A
642	500 mL	N/A	N/A
643	1 L	N/A	N/A
661	10 L	14T8	10/2016

<b>Specifications</b>	<b>Criteria</b>	<b>Lot Data</b>	<b>Results</b>
pH	7.2-7.6	7.38	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	13,320	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control 

Tracy Hanson

Date 10/03/2014

10/03/2014