

CERTIFICATE OF ANALYSIS

SynBlock (BB3)			
Catalog#	Size	Lot#	Expiry
512	Bulk	14B40	11/2016
641	100 mL	N/A	N/A
642	500 mL	N/A	N/A
643	1 L	N/A	N/A
661	10 L	14B41	11/2016

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.34	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMH0s	11,800-16,100	13,950	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Hanson
 Tracy Hanson

Date 11/14/2014
 11/14/2014