

CERTIFICATE OF ANALYSIS

SynBlock (BB3)			
Catalog#	Size	Lot#	Expiry
512	Bulk	17L14	07/2019
641	100 mL	NA	NA
642	500 mL	17L15	07/2019
643	1 L	17L16	07/2019
661	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.33	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	14,550	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy
 Tracy Murphy

Date 07/10/2017
 07/10/2017