

CERTIFICATE OF ANALYSIS

SynBlock (BB3)			
Catalog#	Size	Lot#	Expiry
512	Bulk	18A18	04/2020
641	100 mL	NA	NA
642	500 mL	NA	NA
643	1 L	NA	NA
661	10 L	18A19	04/2020

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.36	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	14,520	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy
 Tracy Murphy

Date 4/10/2018
 4/10/2018