

CERTIFICATE OF ANALYSIS

SynBlock (BB3)			
Catalog#	Size	Lot#	Expiry
512	Bulk	20S19	09/2022
641	100 mL	NA	NA
642	500 mL	NA	NA
643	1 L	NA	NA
661	10 L	20S20	09/2022

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.29	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity μ MHOs	11,800-16,100	14,370	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy
 Tracy Murphy

Date 9/16/2020
 9/16/2020