

CERTIFICATE OF ANALYSIS

SynBlock (BB3)			
Catalog#	Size	Lot#	Expiry
512	Bulk	19F50	02/2021
641	100 mL	NA	NA
642	500 mL	NA	NA
643	1 L	19F51	02/2021
661	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.39	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	14,390	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

Tracy Murphy
 Tracy Murphy

Date

2/26/2019
 2/26/2019