

CERTIFICATE OF ANALYSIS

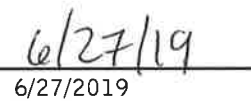
SynBlock (BB3)			
Catalog#	Size	Lot#	Expiry
512	Bulk	18G19	08/2020
641	100 mL	18G20 19N55	08/2020 08/2020
642	500 mL	NA	NA
643	1 L	18G21	08/2020
661	10 L	18G22	08/2020

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.28	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	14,240	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


 Tracy Murphy

Date


 6/27/2019