

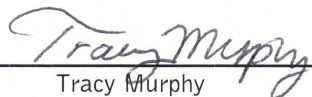
# CERTIFICATE OF ANALYSIS

## SynBlock (BB3)

Catalog#	Size	Lot#	Expiry
512	Bulk	20G33	08/2022
641	100 mL	NA	Pass
642	500 mL	NA	Pass
643	1 L	NA	Pass
661	10 L	20G34	08/2022

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.30	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	14,250	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

  
Tracy Murphy

Date

8/31/2020  
8/31/2020