

CERTIFICATE OF ANALYSIS

SynBlock (BB3)			
Catalog#	Size	Lot#	Expiry
512	Bulk	20N28	06/2022
641	100 mL	NA	NA
642	500 mL	NA	NA
643	1 L	NA	NA
661	10 L	20N29	06/2022

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.27	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	14,620	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control *Sally Hed Dahlquist*
 Sally Hed Dahlquist

Date *6/17/2020*
 6/17/2020