

CERTIFICATE OF ANALYSIS

SynBlock (BB3)			
Catalog#	Size	Lot#	Expiry
512	Bulk	16G57	08/2018
641	100 mL	NA	NA
642	500 mL	NA	NA
643	1 L	16G58	08/2018
661	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.46	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 Tracy Murphy
 Tracy Murphy

Date 08/26/2016
 08/26/2016