

# CERTIFICATE OF ANALYSIS

## SynBlock (BB3)

Catalog#	Size	Lot#	Expiry
512	Bulk	16L5	07/2018
641	100 mL	NA	NA
642	500 mL	NA	NA
643	1 L	NA	NA
661	10 L	16L6	07/2018

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.42	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	14,900	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

*Tracy Murphy*  
 Tracy Murphy

Date

07/06/2016  
 07/06/2016