

# CERTIFICATE OF ANALYSIS

## SynBlock (BB3)

Catalog#	Size	Lot#	Expiry
512	Bulk	15N58	06/2017
641	100 mL	15B6 15D29	06/2017 06/2017
642	500 mL	15N59	06/2017
643	1 L	15N60	06/2017
661	10 L	N/A	N/A

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.32	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	11,800-16,100	14,020	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

*Tracy Murphy*

Tracy Murphy

Date

12/15/2015

12/15/2015