

CERTIFICATE OF ANALYSIS

SynBlock (BB3)			
Catalog#	Size	Lot#	Expiry
512	Bulk	15B19	11/2017
641	100 mL	NA	NA
642	500 mL	NA	NA
643	1 L	NA	NA
661	10 L	15B20	11/2017

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.28	Approved
Appearance	Clear liquid	Conforms	Approved
Conductivity uMHOs	11,800-16,100	13,460	Approved
Storage	Refrigerated	Refrigerated	Approved
Documentation	Approved	Approved	Approved

Quality Control


 Tracy Murphy

Date

11/13/2015
 11/13/2015