

CERTIFICATE OF ANALYSIS

Alternative Block			
Catalog#	Size	Lot#	Expiry
5150	Bulk	15G37	08/2017
6299	100 mL	15G38	08/2017
6300	500 mL	N/A	N/A
6301	1 L	15G39	08/2017
6302	10 L	N/A	N/A

Specifications	Criteria	Lot Data	Results
pH	7.1-7.5	7.46	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity μ MHOs	Report measurement	14,630	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Hanson
 Tracy Hanson

Date 08/18/2015
 08/18/2015