

CERTIFICATE OF ANALYSIS

Alternate Block™			
Catalog#	Size	Lot#	Expiry at Room Temperature
5150	Bulk	19Y40	05/2021
6299	100 mL	19Y41	05/2021
6301	1 L	19Y42	05/2021
6302	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.1-7.6	7.41	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity μ MHOs	13,000-20,000	15,340	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

This is an assurance that the specified lot conforms to all requirements of the product specifications.
 This information is believed to be accurate and is offered in good faith.
 Product is for research use only. Not for use in diagnostic procedures.

Quality Assurance: 
 Tracy Murphy

Date: 5/16/2019
 5/16/2019