



# CERTIFICATE OF ANALYSIS

Protein-Free Sample Diluent			
Catalog#	Size	Lot#	Expiry at 2-8°C
5178	Bulk	20A14	04/2022
6702	100 mL	20A15	04/2022
6703	500 mL	20A16	04/2022
6704	1 L	20A17	04/2022
6707	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.1-7.6	7.34	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity $\mu$ MHOs	13,000-20,000	16,820	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.

Quality Assurance *Sally Hed Dahlquist*  
Sally Hed Dahlquist

Date 4/20/2020  
4/20/2020

9401 James Avenue S., #155  
Bloomington, MN 55431 USA  
TOLL-FREE 1-800-829-3194  
LOCAL 952-888-8788