



# CERTIFICATE OF ANALYSIS

Protein-Free Sample Diluent			
Catalog#	Size	Lot#	Expiry at 2-8°C
5178	Bulk	18B28	11/2020
6702	100 mL	18B29	11/2020
6703	500 mL	18B30	11/2020
6704	1 L	18B31	11/2020
6707	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
<b>pH</b>	7.1-7.6	7.47	Pass
<b>Appearance</b>	Clear liquid	Conforms	Pass
<b>Conductivity <math>\mu</math>MHOs</b>	13,000-20,000	18,570	Pass
<b>Storage</b>	Refrigerated	Refrigerated	Pass
<b>Documentation</b>	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   
 Kristi Strandberg

Date 11/13/18  
 11/13/2018

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

IMMUNOCHEMISTRY.COM