

# CERTIFICATE OF ANALYSIS

## General Assay Diluent (AD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
518	Bulk	16S11	09/2018
620	100 mL	NA	NA
621	500 mL	NA	NA
622	1 L	NA	NA
671	10 L	16S12	09/2018

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.43	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMHOs	40,000-60,000	43,600	Pass
A <sub>280</sub> of 1:20 dilution	0.225-0.375	0.291	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

*Tracy Murphy*  
 Tracy Murphy

Date

09/12/2016  
 09/12/2016