

# CERTIFICATE OF ANALYSIS

## General Assay Diluent (AD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
518	Bulk	19R42	03/2021
620	100 mL	19R43	03/2021
621	500 mL	19R44	03/2021
622	1 L	19R45 19Y9	03/2021 03/2021
671	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.26	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMHOs	40,000-60,000	41,800	Pass
A <sub>280</sub> of 1:20 dilution	0.225-0.375	0.260	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control



Kristi Strandberg

Date

05/03/19

5/3/2019