

CERTIFICATE OF ANALYSIS

General Assay Diluent (AD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
518	Bulk	19G55	08/2021
620	100 mL	NA	NA
621	500 mL	NA	NA
622	1 L	19G56	08/2021
671	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.41	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMHOs	40,000-60,000	42,500	Pass
A ₂₈₀ of 1:20 dilution	0.225-0.375	0.265	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

Tracy Murphy
Tracy Murphy

Date

8/22/19
8/22/2019