

CERTIFICATE OF ANALYSIS

General Assay Diluent (AD1)


Catalog#	Size	Lot#	Expiry at 2-8°C
518	Bulk	18A23	04/2020
620	100 mL	NA	NA
621	500 mL	NA	NA
622	1 L	18A24	04/2020
671	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.39	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMHOs	40,000-60,000	40,010	Pass
A ₂₈₀ of 1:20 dilution	0.225-0.375	0.271	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


 Tracy Murphy

Date


 4/16/2018