

CERTIFICATE OF ANALYSIS

General Assay Diluent (AD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
518	Bulk	15T8	10/2017
620	100 mL	16Y6	10/2017
621	500 mL	15T9	10/2017
622	1 L	15T10	10/2017
671	10 L	N/A	N/A

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.30	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMH0s	40,000-60,000	59,700	Pass
A ₂₈₀ of 1:20 dilution	0.225-0.375	0.246	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

Tracy Murphy

Tracy Murphy

Date

05/06/2016

05/06/2016