

CERTIFICATE OF ANALYSIS

General Serum Diluent (SD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
526	Bulk	18Y23	05/2020
647	100 mL	NA	NA
648	500 mL	NA	NA
649	1 L	18Y24	05/2020
675	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.27	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMH0s	14,000-21,000	20,100	Pass
A ₂₈₀ of 1:20 dilution	0.48-0.72	0.685	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

Tracy Murphy
 Tracy Murphy

Date

05/21/2018
 05/21/2018