

CERTIFICATE OF ANALYSIS

General Serum Diluent (SD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
526	Bulk	17D24	12/2019
647	100 mL	NA	NA
648	500 mL	NA	NA
649	1 L	NA	NA
675	10 L	17D25	12/2019

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.31	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMHOs	14,000-21,000	17,860	Pass
A ₂₈₀ of 1:20 dilution	0.48-0.72	0.623	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


 Tracy Murphy

Date

12/09/2017
 12/09/2017