

CERTIFICATE OF ANALYSIS

General Serum Diluent (SD1)

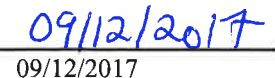
Catalog#	Size	Lot#	Expiry at 2-8°C
526	Bulk	17S10	09/2019
647	100 mL	NA	NA
648	500 mL	NA	NA
649	1 L	17S11	09/2019
675	10 L	NA	NA

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	17,910	Pass
A ₂₈₀ of 1:20 dilution	0.48-0.72	0.628	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


 Tracy Murphy

Date


 09/12/2017