

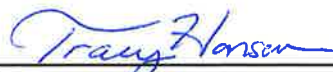
# CERTIFICATE OF ANALYSIS

## General Serum Diluent (SD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
526	Bulk	15R57	03/2017
647	100 mL	N/A	N/A
648	500 mL	N/A	N/A
649	1 L	15R58	03/2017
675	10 L	N/A	N/A

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.36	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMH0s	14,000-21,000	16,260	Pass
A <sub>280</sub> of 1:20 dilution	0.48-0.72	0.638	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control



Tracy Hanson

Date

03/18/2015

03/18/2015