

# CERTIFICATE OF ANALYSIS

## General Serum Diluent (SD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
526	Bulk	14S1	09/2016
647	100 mL	N/A	N/A
648	500 mL	N/A	N/A
649	1 L	14S2	09/2016
675	10 L	N/A	N/A

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.35	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMH0s	14,000-21,000	17,120	Pass
A <sub>280</sub> of 1:20 dilution	0.48-0.72	0.635	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Hanson

Tracy Hanson

Date 09/05/2014

09/05/2014