

# CERTIFICATE OF ANALYSIS

## General Serum Diluent (SD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
526	Bulk	16S26	09/2018
647	100 mL	16S27	09/2018
648	500 mL	16S28	09/2018
649	1 L	16S29	09/2018
675	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.36	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMHOs	14,000-21,000	17,920	Pass
A <sub>280</sub> of 1:20 dilution	0.48-0.72	0.642	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

  
 Tracy Murphy

Date

09/23/2016  
 09/23/2016