

# CERTIFICATE OF ANALYSIS

## General Serum Diluent (SD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
526	Bulk	17L31	07/2019
647	100 mL	NA	NA
648	500 mL	NA	NA
649	1 L	17L32	07/2019
675	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.37	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMH0s	14,000-21,000	18,060	Pass
A <sub>280</sub> of 1:20 dilution	0.48-0.72	0.620	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy  
 Tracy Murphy

Date 07/25/2017  
 07/25/2017