

# CERTIFICATE OF ANALYSIS

## General Serum Diluent (SD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
526	Bulk	20Y6	05/2022
647	100 mL	NA	NA
648	500 mL	NA	NA
649	1 L	NA	NA
675	10 L	20Y7	05/2022

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.45	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMHOs	14,000-21,000	18,090	Pass
A <sub>280</sub> of 1:20 dilution	0.48-0.72	0.596	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

  
 Sally Hed Dahlquist

Date

  
 5/7/2020