


CERTIFICATE OF ANALYSIS

General Serum Diluent (SD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
526	Bulk	16A10	04/2018
647	100 mL	16L27	04/2018
648	500 mL	16A11	04/2018
649	1 L	16A12	04/2018
675	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.34	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMH0s	14,000-21,000	17,960	Pass
A ₂₈₀ of 1:20 dilution	0.48-0.72	0.660	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


 Tracy Murphy

Date

07/19/2016
 07/19/2016