

CERTIFICATE OF ANALYSIS

General Serum Diluent (SD1)

Catalog#	Size	Lot#	Expiry at 2-8°C
526	Bulk	14L33	07/2016
647	100 mL	NA	NA
648	500 mL	NA	NA
649	1 L	14L34	07/2016
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	16,260	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 Hanson

Date

07/21/2014
07/21/2014