

CERTIFICATE OF ANALYSIS

General Serum Diluent (SD1)

Catalog #	Size	Lot #	Expiry at 2-8°C
526	Bulk	14B54	11/2016
647	100 mL	N/A	N/A
648	500 mL	N/A	N/A
649	1 L	14B55	11/2016
675	10 L	N/A	N/A

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.35	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMH0s	14,000-21,000	16,370	Pass
A ₂₈₀ of 1:20 dilution	0.48-0.72	0.628	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control



Tracy Hanson

Date

11/19/2014

11/19/2014