

CERTIFICATE OF ANALYSIS

General Serum Diluent (SD1)			
Catalog#	Size	Lot#	Expiry at 2-8°C
526	Bulk	15Y45	05/2017
647	100 mL	NA	NA
648	500 mL	NA	NA
649	1 L	15Y46	05/2017
675	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.42	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMHOs	14,000-21,000	16,440	Pass
A ₂₈₀ of 1:20 dilution	0.48-0.72	0.659	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Hanson

Date 06/01/2015

Tracy Hanson

06/01/2015