

CERTIFICATE OF ANALYSIS

Plasma Sample Diluent (SD2)

Catalog#	Size	Lot#	Expiry at 2-8°C
5123	Bulk	19A39	04/2021
694	100 mL	19A40	04/2021
695	500 mL	19A41	04/2021
696	1 L	19A42	04/2021
697	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.40	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity μ MHOs	30,000-55,000	31,900	Pass
A ₂₈₀ of 1:20 dilution	0.28-0.42	0.344	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

Tracy Murphy
 Tracy Murphy

Date

4/29/19
 4/29/2019