

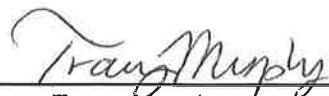
CERTIFICATE OF ANALYSIS

Plasma Sample Diluent (SD2)

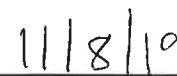
Catalog#	Size	Lot#	Expiry at 2-8°C
5123	Bulk	19B10	11/2021
694	100 mL	19B11	11/2021
695	500 mL	19B12	11/2021
696	1 L	19B13	11/2021
697	10 L	NA	NA

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	32,500	Pass
A ₂₈₀ of 1:20 dilution	0.28-0.42	0.342	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


Tracy Murphy

Date


11/8/2019