

# CERTIFICATE OF ANALYSIS

## IgM-Reducing Assay Diluent (AD2)

Catalog#	Size	Lot#	Expiry at 2-8°C
519	Bulk	19B26	05/2021
623	100 mL	20A8	05/2021
624	500 mL	19B27	05/2021
625	1 L	19B28	05/2021
672	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	6.4-6.6	6.48	Pass
Appearance	Clear to cloudy, slightly white liquid	Conforms	Pass
Conductivity uMHOs	56,000-84,000	69,500	Pass
A <sub>280</sub> of 1:20 dilution	0.12-0.18	0.164	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

*Sally Dahlquist*  
Sally Dahlquist

Date

*4/17/2020*  
04/17/2020