

CERTIFICATE OF ANALYSIS

IgM-Reducing Assay Diluent (AD2)

Catalog#	Size	Lot#	Expiry at 2-8°C
519	Bulk	19G43	02/2021
623	100 mL	NA	NA
624	500 mL	NA	NA
625	1 L	19G44	02/2021
672	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	6.4-6.6	6.45	Pass
Appearance	Clear to cloudy, slightly white liquid	Conforms	Pass
Conductivity uMH0s	56,000-84,000	61,200	Pass
A ₂₈₀ of 1:20 dilution	0.12-0.18	0.172	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy
 Tracy Murphy

Date 8/21/19
 8/21/2019