

CERTIFICATE OF ANALYSIS

IgM-Reducing Assay Diluent (AD2)

Catalog#	Size	Lot#	Expiry at 2-8°C
519	Bulk	20N22	12/2021
623	100 mL	20N23	12/2021
624	500 mL	20N24	12/2021
625	1 L	20N25	12/2021
672	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	6.4-6.6	6.47	Pass
Appearance	Clear to cloudy, slightly white liquid	Conforms	Pass
Conductivity uMHOs	56,000-84,000	67,500	Pass
A ₂₈₀ of 1:20 dilution	0.12-0.18	0.156	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


Sally Hed Dahlquist

Date


6/12/2020