

CERTIFICATE OF ANALYSIS

IgM-Reducing Assay Diluent (AD2)

Catalog#	Size	Lot#	Expiry at 2-8°C
519	Bulk	17A16	10/2018
623	100 mL	17A17	10/2018
624	500 mL	17A18	10/2018
625	1 L	17A19	10/2018
672	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	6.4-6.6	6.51	Pass
Appearance	Clear to cloudy, slightly white liquid	Conforms	Pass
Conductivity uMHOs	56,000-84,000	61,300	Pass
A ₂₈₀ of 1:20 dilution	0.12-0.18	0.151	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control 
 Kristi Strandberg

Date 04/11/17
 04/11/2017