

CERTIFICATE OF ANALYSIS

IgM-Reducing Assay Diluent (AD2)

Catalog#	Size	Lot#	Expiry at 2-8°C
519	Bulk	15T11	10/2017
623	100 mL	15T12	10/2017
624	500 mL	15T13	10/2017
625	1 L	15T14	10/2017
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 uMH0s	56,000-84,000	74,800	Pass
A ₂₈₀ of 1:20 dilution	0.12-0.18	0.164	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

Tracy Murphy
 Tracy Murphy

Date

10/12/2015
 10/12/2015