

# CERTIFICATE OF ANALYSIS

## IgM-Reducing Assay Diluent (AD2)

Catalog#	Size	Lot#	Expiry at 2-8°C
519	Bulk	14B62	11/2016
623	100 mL	14B63 15L36	11/2016 11/2016
624	500 mL	14B64	11/2016
625	1 L	14B65	11/2016
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	59,700	Pass
A <sub>280</sub> of 1:20 dilution	0.12-0.18	0.136	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

  
 Tracy Hanson

Date

07/17/2015  
 07/17/2015