

CERTIFICATE OF ANALYSIS

IgM-Reducing Assay Diluent (AD2)

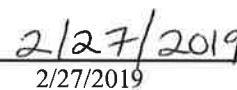
| Catalog# | Size | Lot# | Expiry at 2-8°C |
|----------|--------|-------|-----------------|
| 519 | Bulk | 19F53 | 08/2020 |
| 623 | 100 mL | NA | NA |
| 624 | 500 mL | 19F54 | 08/2020 |
| 625 | 1 L | NA | NA |
| 672 | 10 L | NA | NA |

| Specifications | Criteria | Lot Data | Results |
|-----------------------------------|--|--------------|---------|
| pH | 6.4-6.6 | 6.48 | 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 | 0.144 | Pass |
| Storage | Refrigerated | Refrigerated | Pass |
| Documentation | Approved | Approved | Pass |

Quality Control


Tracy Murphy

Date


2/27/2019