

CERTIFICATE OF ANALYSIS

General Serum Diluent (SD1)

| Catalog# | Size | Lot# | Expiry at 2-8°C |
|----------|--------|-------|-----------------|
| 526 | Bulk | 17N9 | 06/2019 |
| 647 | 100 mL | NA | NA |
| 648 | 500 mL | NA | NA |
| 649 | 1 L | 17N10 | 06/2019 |
| 675 | 10 L | NA | NA |

| Specifications | Criteria | Lot Data | Results |
|-----------------------------------|---------------------------------|--------------|---------|
| pH | 7.2-7.6 | 7.24 | Pass |
| Appearance | Clear to slightly yellow liquid | Conforms | Pass |
| Conductivity uMH0s | 14,000-21,000 | 18,120 | Pass |
| A ₂₈₀ of 1:20 dilution | 0.48-0.72 | 0.612 | Pass |
| Storage | Refrigerated | Refrigerated | Pass |
| Documentation | Approved | Approved | Pass |

Quality Control

Tracy Murphy
 Tracy Murphy

Date

06/08/2017
 06/08/2017