

CERTIFICATE OF ANALYSIS

| General Serum Diluent (SD1) | | | |
|------------------------------------|-------------|-------------|------------------------|
| Catalog# | Size | Lot# | Expiry at 2-8°C |
| 526 | Bulk | 15R86 | 03/2017 |
| 647 | 100 mL | 15R87 | 03/2017 |
| 648 | 500 mL | 15R88 | 03/2017 |
| 649 | 1 L | N/A | N/A |
| 675 | 10 L | N/A | N/A |

| Specifications | Criteria | Lot Data | Results |
|-----------------------------------|---------------------------------|-----------------|----------------|
| pH | 7.2-7.6 | 7.28 | Pass |
| Appearance | Clear to slightly yellow liquid | Conforms | Pass |
| Conductivity uMHOs | 14,000-21,000 | 16,450 | Pass |
| A ₂₈₀ of 1:20 dilution | 0.48-0.72 | 0.629 | Pass |
| Storage | Refrigerated | Refrigerated | Pass |
| Documentation | Approved | Approved | Pass |

Quality Control



Tracy Hanson

Date

03/26/2015

03/26/2015