

CERTIFICATE OF ANALYSIS

| Stop Solution for TMB Substrates (STOPT) | | | |
|--|--------|-------|------------|
| Catalog# | Size | Lot# | Expiry |
| 3331 | Bulk | 17J23 | 07/28/2019 |
| 6282 | 100 mL | NA | NA |
| 6343 | 1 L | 17J24 | 07/28/2019 |

| Specifications | Criteria | Lot Data | Results |
|----------------|---------------------------|------------------|---------|
| Appearance | Clear, Colorless Solution | Conforms | Pass |
| pH | ≤ 2.00 | 1.20 | Pass |
| Storage | Room Temperature (25°C) | Room Temperature | Pass |

This is an assurance that the specified lot conforms to all requirements of the product specifications.
 This information is believed to be accurate and is offered in good faith.
 Product is for research use only. Not for use in diagnostic procedures.

Quality Assurance: Tracy Murphy
 Tracy Murphy

Date: 01/17/2017
 01/17/2017