



CERTIFICATE OF ANALYSIS

Stop Solution for TMB Substrates (STOPT)			
Catalog#	Size	Lot#	Expiry at 25°C
3331	Bulk	19R49	12/04/2021
6282	100 mL	19A24	12/04/2021
6343	1 L	19R50	12/04/2021

Specifications	Criteria	Lot Data	Results
pH	≤ 2.00	Conforms	Pass
Appearance	Clear, colorless	1.19	Pass
Storage	Room temperature (25°C)	Room Temperature (25°C)	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.

Quality Assurance Tracy Murphy
 Tracy Murphy

Date 4/15/2019
 4/15/2019