



CERTIFICATE OF ANALYSIS

Stop Solution for TMB Substrates (STOPT)			
Catalog#	Size	Lot#	Expiry at 25°C
3331	Bulk	19N63	01/18/2022
6282	100 mL	NA	NA
6343	1 L	19N64	01/18/2022

Specifications	Criteria	Lot Data	Results
pH	≤ 2.00	1.21	Pass
Appearance	Clear, colorless	Conforms	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 7/1/19
 7/01/2019