



# CERTIFICATE OF ANALYSIS

Stop Solution for TMB Substrates (STOPT)			
Catalog#	Size	Lot#	Expiry at 25°C
3331	Bulk	19Y64	01/18/2022
6282	100 mL	NA	NA
6343	1 L	19Y65	01/18/2022

Specifications	Criteria	Lot Data	Results
<b>pH</b>	≤ 2.00	1.21	Pass
<b>Appearance</b>	Clear, colorless	Conforms	Pass
<b>Storage</b>	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 5/30/19  
5/30/2019