



CERTIFICATE OF ANALYSIS

Stop Solution for TMB Substrates (STOPT)			
Catalog#	Size	Lot#	Expiry at 25°C
3331	Bulk	21J5	12/10/2023
6282	100 mL	NA	NA
6343	1 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	≤ 2.00	1.25	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 _____ *1/11/2021*
 1/11/2021