

# CERTIFICATE OF ANALYSIS

Stop Solution for TMB Substrates (STOPT)			
Catalog#	Size	Lot#	Expiry
3331	Bulk	17J45	01/20/2020
6282	100 mL	NA	NA
6343	1 L	NA	NA

Specifications	Criteria	Lot Data	Results
Appearance	Clear, Colorless Solution	Conforms	Pass
pH	≤ 2.00	1.22	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: 01/27/2017  
 01/27/2017