

CERTIFICATE OF ANALYSIS

Stop Solution for TMB Substrates (STOPT)			
Catalog#	Size	Lot#	Expiry
3331	Bulk	16F11	06/18/2018
6282	100 mL	NA	NA
6343	1 L	NA	NA

Specifications	Criteria	Lot Data	Results
Appearance	Clear, Colorless Solution	Conforms	Pass
pH	≤ 2.00	1.19	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.

Signature: 
 Quality Assurance

Date: 02/05/2016
 02/05/2016