

# CERTIFICATE OF ANALYSIS

TMB One Component HRP Microwell Substrate (SUB1)			
Catalog#	Size	Lot#	Expiry
3325	Bulk	13B39	08/09/2016
6276	100 mL	13B40 14A28	08/09/2016 08/09/2016

Specifications	Criteria	Lot Data	Results
pH	3.55 ± 0.20	3.62	Pass
Absorbance at 285 nm	0.450 ± 0.100	0.439	Pass
Absorbance at 650 nm	≤ 0.020	0.000	Pass
Appearance	Colorless to very light yellow solution	Conforms	Pass
Storage	Refrigerated	Refrigerated	Pass

Quality Control Tracy Hanson

Tracy Hanson

Date 04/21/2014

04/21/2014