


CERTIFICATE OF ANALYSIS

TMB Super Sensitive HRP Microwell Substrate (SUB2)			
Catalog#	Size	Lot#	Expiry
3324	Bulk	14A34	08/03/2016
6275	100 mL	NA	NA
6328	500 mL	NA	NA
6329	1 L	NA	NA
6330	10 L	14A35	08/03/2016

Specifications	Criteria	Lot Data	Results
pH	3.70 ± 0.20	3.75	Pass
Absorbance at 285 nm	0.600 ± 0.150	0.524	Pass
Absorbance at 650 nm	≤ 0.050	0.002	Pass
Appearance	Colorless to very light yellow solution	Conforms	Pass
Storage	Refrigerated	Refrigerated	Pass

Quality Control



Tracy Hanson

Date

04/21/2014

04/21/2014