

CERTIFICATE OF ANALYSIS

1X HRP Conjugate Stabilizer Diluent			
Catalog#	Size	Lot#	Expiry at 2-8°C
5170	Bulk	17Y35	05/2019
6350	100 mL	N/A	N/A
6351	1 L	N/A	N/A
6352	10 L	17Y36	05/2019

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.24	Pass
Appearance	Clear to slightly cloudy, yellow tinted liquid	Conforms	Pass
Conductivity	Report Results	16,460 μ MHOs	Pass
Spec A ₂₈₀ (1:20)	Report Results	0.902	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control



Kristi Strandberg

Date

05/25/17

05/25/2017