

CERTIFICATE OF ANALYSIS

1X HRP Conjugate Stabilizer Diluent			
Catalog#	Size	Lot#	Expiry at 2-8°C
5170	Bulk	19N33	06/2021
6350	100 mL	NA	NA
6351	1 L	19N34	06/2021
6352	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.35	Pass
Appearance	Clear to slightly cloudy, yellow tinted liquid	Conforms	Pass
Conductivity	Report Results	16,060	Pass
Spec A ₂₈₀ (1:20)	Report Results	0.870	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

Tracy Murphy
 Tracy Murphy

Date

6/14/19
 6/14/2019