

CERTIFICATE OF ANALYSIS

1X HRP Conjugate Stabilizer Diluent

Catalog#	Size	Lot#	Expiry at 2-8°C
5170	Bulk	15S14	09/2017
6350	100 mL	15S15	09/2017
6351	1 L	15S16	09/2017
6352	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.40	Pass
Appearance	Clear to slightly cloudy, yellow tinted liquid	Conforms	Pass
Conductivity	Report Results	15,050	Pass
Spec A ₂₈₀ (1:20)	Report Results	0.955	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

Tracy Hanson
 Tracy Hanson

Date

09/09/2015
 09/09/2015