

CERTIFICATE OF ANALYSIS

1X Neptune HRP Conjugate Stabilizer Diluent			
Catalog#	Size	Lot#	Expiry at 2-8°C
5169	Bulk	19L10	07/2021
6347	100 mL	19L11	07/2021
6348	1 L	NA	NA
6349	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.39	Pass
Appearance	Clear to slightly cloudy, yellow tinted liquid	Conforms	Pass
Conductivity	Report Results	16,400 μ MHOs	Pass
Spec A ₂₈₀ (1:20)	Report Results	0.589	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


 Kristi Strandberg

Date

07/03/19

07/03/2019