



CERTIFICATE OF ANALYSIS

Alkaline Phosphatase Conjugate Stabilizer, 1X			
Catalog#	Size	Lot#	Expiry at 2-8°C
3323	Bulk	20J2	12/03/2021
6270	25 mL	NA	NA
6271	100 mL	NA	NA
6273	1 L	NA	NA
6274	10 L	20J3	12/03/2021

Specifications	Criteria	Lot Data	Results
pH	6.0 ± 0.50	6.0	Pass
Accelerated Stability Test: AP antibody conjugate is diluted to working concentration, stored at 37°C for 1 week, and then tested for retained activity.	≥85% (relative to 4°C control)	Conforms	Pass
Appearance	Clear, light yellow liquid	Conforms	Pass
Storage	Refrigerated	Refrigerated	Pass

This is an assurance that the specified lot conforms to all requirements of the product specifications. This information is believed to be accurate and is offered in good faith.

Quality Assurance Tracy Murphy
Tracy Murphy

Date 1/3/2020
1/3/2020