

CERTIFICATE OF ANALYSIS

Alkaline Phosphatase Conjugate Stabilizer, 1X			
Catalog#	Size	Lot#	Expiry
3323	Bulk	NA	NA
6270	25 mL	NA	NA
6271	100 mL	15R35	07/29/2016
6273	1 L	14N57	07/29/2016
6274	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	6.0 ± 0.50	Conforms	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.

Signature: Tracy Hanson

Tracy Hanson

Date: 03/18/2015

03/18/2015