

CERTIFICATE OF ANALYSIS

5X Antibody Coating Buffer (CB1)			
Catalog#	Size	Lot#	Expiry at room temperature
524	Bulk	14D26	12/2016
6245	25 mL	N/A	N/A
644	100 mL	N/A	N/A
645	500 mL	15A25	12/2016
646	1 L	14D27	12/2016
658	10 L	N/A	N/A

Specifications	Criteria	Lot Data	Results
pH at 5X	7.8-8.2	8.10	Pass
pH at 1X	8.2-8.6	8.20	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs at 5X	48,000-65,000	64,900	Pass
Storage	Refrigerated or room temperature	Room Temperature	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Hanson

Date 04/10/2014

Tracy Hanson

04/10/2014