

CERTIFICATE OF ANALYSIS

5X Antibody Coating Buffer (CB1)			
Catalog#	Size	Lot#	Expiry at room temperature
524	Bulk	16G46	08/2018
6245	25 mL	NA	NA
644	100 mL	16B44	08/2018
645	500 mL	16S18	08/2018
646	1 L	16G47	08/2018
658	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH at 5X	7.8-8.2	8.10	Pass
pH at 1X	8.2-8.6	8.24	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs at 5X	48,000-65,000	49,200	Pass
Storage	Refrigerated or room temperature	Room temperature	Pass
Documentation	Approved	Approved	Pass

Quality Control 
Tracy Murphy

Date 11/29/2016
11/29/2016