

CERTIFICATE OF ANALYSIS

5X Antibody Coating Buffer (CB1)			
Catalog#	Size	Lot#	Expiry at room temperature
524	Bulk	16N28	06/2018
6245	25 mL	NA	NA
644	100 mL	NA	NA
645	500 mL	16N59	06/2018
646	1 L	16N29	06/2018
658	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH at 5X	7.8-8.2	8.14	Pass
pH at 1X	8.2-8.6	8.39	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMH0s at 5X	48,000-65,000	50,800	Pass
Storage	Refrigerated or room temperature	Room temperature	Pass
Documentation	Approved	Approved	Pass

Quality Control


Tracy Murphy

Date

06/23/2016
06/23/2016