

# CERTIFICATE OF ANALYSIS

## 5X Antibody Coating Buffer (CB1)

Catalog#	Size	Lot#	Expiry at room temperature
524	Bulk	15D1	12/2017
6245	25 mL	NA	NA
644	100 mL	NA	NA
645	500 mL	15D2	12/2017
646	1 L	NA	NA
658	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH at 5X	7.8-8.2	8.18	Pass
pH at 1X	8.2-8.6	8.34	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs at 5X	48,000-65,000	56,600	Pass
Storage	Refrigerated or room temperature	Room Temperature	Pass
Documentation	Approved	Approved	Pass

Quality Control

*Tracy Murphy*  
 Tracy Murphy

Date

*12/02/2017*  
 12/02/2017