

CERTIFICATE OF ANALYSIS

5X Antigen Coating Buffer (CB2)

Catalog#	Size	Lot#	Expiry at room temperature
5136	Bulk	16G48	08/2018
6246	25 mL	NA	NA
6247	100 mL	16G49	08/2018
6248	500 mL	NA	NA
6249	1 L	16G50	08/2018
6250	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

08/25/2016
08/25/2016