

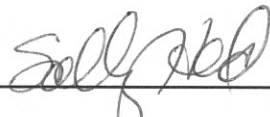
# CERTIFICATE OF ANALYSIS

## Neptune Blocking Buffer (BB2)

Catalog#	Size	Lot#	Expiry
53	Bulk	14T52	10/2016
62	100 mL	14T72	10/2016
63	500 mL	14T53	10/2016
64	1 L	N/A	N/A
660	10 L	14T54	10/2016

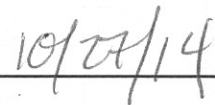
Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.6	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity uMHOs	15,000-19,000	15,610	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control



Sally Hed

Date



10/27/2014