

CERTIFICATE OF ANALYSIS

Neptune Blocking Buffer (BB2)

Catalog#	Size	Lot#	Expiry
53	Bulk	15F23	02/2017
62	100 mL	15F24	02/2017
63	500 mL	15F25	02/2017
64	1 L	15F26	02/2017
660	10 L	N/A	N/A

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.44	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity uMH0s	15,000-19,000	15,230	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Hanson

Tracy Hanson

Date 02/13/2015

02/13/2015