

CERTIFICATE OF ANALYSIS

Neptune Blocking Buffer (BB2)

Catalog#	Size	Lot#	Expiry
53	Bulk	15S10	09/2017
62	100 mL	16N76	NA
63	500 mL	16J42 16N75	09/2017 09/2017
64	1 L	15S11	09/2017
660	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.43	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity uMH0s	15,000-19,000	15,340	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy
 Tracy Murphy

Date 07/20/16
 07/20/2016