

CERTIFICATE OF ANALYSIS

Neptune Blocking Buffer (BB2)

Catalog#	Size	Lot#	Expiry
53	Bulk	17L24	07/2019
62	100 mL	NA	NA
63	500 mL	NA	NA
64	1 L	17L25	07/2019
660	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.50	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity uMHOs	15,000-19,000	16,520	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy
 Tracy Murphy

Date 07/21/2017
 07/21/2017