



CERTIFICATE OF ANALYSIS

Neptune Block			
Catalog#	Size	Lot#	Expiry at 2-8°C
53	Bulk	21J2	01/2023
62	100 mL	NA	NA
63	500 mL	NA	NA
64	1 L	21J3	01/2023
660	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.42	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity μMHOs	15,000-19,000	16,440	Pass
A₂₈₀ of 1:10 dilution	0.34-0.68	0.541	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

This is an assurance that the specified lot conforms to all requirements of the product specifications. This information is believed to be accurate and is offered in good faith.

Quality Assurance Tracy Murphy
 Tracy Murphy

Date 1/6/2021
 1/06/2021