



CERTIFICATE OF ANALYSIS

Neptune Assay Diluent			
Catalog#	Size	Lot#	Expiry at 2-8°C
522	Bulk	20J35	01/2022
626	100 mL	20J36	01/2022
627	500 mL	20J37	01/2022
628	1 L	20J38 20G8	01/2022 01/2022
673	10 L	N/A	N/A

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.43	Pass
Appearance	Light-yellow liquid	Conforms	Pass
Conductivity μMHOs	15,000-45,000	32,900	Pass
A₂₈₀ of 1:20 dilution	0.2-0.5	0.291	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 Sally Hed Dahlquist
Sally Hed Dahlquist

Date 8/14/2020
8/14/2020

9401 James Avenue S., #155
Bloomington, MN 55431 USA
TOLL-FREE 1-800-829-3194
LOCAL 952-888-8788