

CERTIFICATE OF ANALYSIS

Neptune™ Sample Diluent

Catalog#	Size	Lot#	Expiry at 2-8°C
591	Bulk	20S35	09/2022
6124	100 mL	NA	NA
6125	500 mL	NA	NA
6126	1 L	20S36	09/2022
6127	10 L	NA	NA

Specification	Criteria	Lot Data	Result
pH	7.2-7.6	7.41	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMHOs	Report result	21,800	Pass
A ₂₈₀ of 1:20 dilution	Report result	0.153	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.

Signature: Traug Murphy
 Quality Assurance

Date: 9/25/2020
 9/25/2020