

CERTIFICATE OF ANALYSIS

Neptune™ Sample Diluent			
Catalog#	Size	Lot#	Expiry at 2-8°C
591	Bulk	19L57	07/2021
6124	100 mL	19L58	07/2021
6125	500 mL	19L59	07/2021
6126	1 L	19L60	07/2021
6127	10 L	NA	NA

Specification	Criteria	Lot Data	Result
pH	7.2-7.6	7.40	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity uMH0s	Report result	20,400	Pass
A ₂₈₀ of 1:20 dilution	Report result	0.159	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: Tracy Murphy
 Tracy Murphy

Date: 7/26/19
 7/26/2019