

CERTIFICATE OF ANALYSIS

Neptune Assay Diluent (AD3)

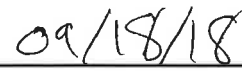
Catalog#	Size	Lot#	Expiry at 2-8°C
522	Bulk	18S26	09/2020
626	100 mL	18S27	09/2020
627	500 mL	NA	NA
628	1 L	18S28	09/2020
673	10 L	NA	NA

Specification	Criteria	Lot Data	Result
pH	7.2-7.6	7.45	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity μ MHO	20,000-50,000	30,600	Pass
A ₂₈₀ of 1:20 dilution	0.2-0.5	0.336	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


 Kristi Standberg

Date


 9/18/2018