

CERTIFICATE OF ANALYSIS

Neptune Assay Diluent (AD3)

Catalog#	Size	Lot#	Expiry at 2-8°C
522	Bulk	15N34	06/2017
626	100 mL	15N35 16R16	06/2017 06/2017
627	500 mL	15N36	06/2017
628	1 L	15N37	06/2017
673	10 L	NA	NA

Specification	Criteria	Lot Data	Result
pH	7.2-7.6	7.26	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity μ MHO	20,000-50,000	43,600	Pass
A ₂₈₀ of 1:20 dilution	0.2-0.5	0.323	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


Tracy Murphy

Date

03/09/2016
03/09/2016