



9401 James Avenue S., Suite #155, Bloomington, MN 55431 USA
952-888-8788 fax 952-888-8988

Document: F11-522-4-C
Effective: 11/13/13
Supersedes: F11-522-4-B
Page 1 of 1

CERTIFICATE OF ANALYSIS

Neptune Assay Diluent (AD3)

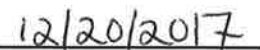
Catalog#	Size	Lot#	Expiry at 2-8°C
522	Bulk	16B6	11/2018
626	100 mL	16B7	11/2018
627	500 mL	16B8 17D52	11/2018 11/2018
628	1 L	16B9	11/2018
673	10 L	NA	NA

Specification	Criteria	Lot Data	Result
pH	7.2-7.6	7.48	Pass
Appearance	Light yellow liquid	Conforms	Pass
Conductivity μ MHO	20,000-50,000	35,700	Pass
A ₂₈₀ of 1:20 dilution	0.2-0.5	0.327	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


Tracy Murphy

Date


12/20/2017