



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

Document: F11-52-4-A
Effective: 10/31/12
Supersedes: None
Page 1 of 1

CERTIFICATE OF ANALYSIS


5X HRP Conjugate Stock Stabilizer Diluent (5X CS6)			
Catalog#	Size	Lot#	Expiry at 2-8°C
52	Bulk	18B46	11/2020
6173	25 mL	19Y18	11/2020
667	100 mL	18B47	11/2020
668	500 mL	18B48	11/2020
669	1 L	18B49	11/2020
670	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH at 5X	7.2-7.6	7.41	Pass
Appearance	Clear to slightly cloudy, yellow tinted liquid	Conforms	Pass
Conductivity uMHOs	60,000-80,000 uMHO	63,400	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


Tracy Murphy

Date


05/07/2019