

CERTIFICATE OF ANALYSIS

5X Antigen-Down HRP Conjugate Stabilizer Diluent (5X CS2)			
Catalog#	Size	Lot#	Expiry at 2-8°C
586	Bulk	16N72	07/2018
6169	25 mL	NA	NA
6102	100 mL	NA	NA
6103	500 mL	16N73	07/2018
6104	1 L	16G37 16T5	07/2018 07/2018
6105	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH at 5X	7.2-7.6	7.4	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
 Tracy Murphy

Date

10/04/2016
 10/04/2016