

CERTIFICATE OF ANALYSIS

5X Antigen-Down HRP Conjugate Stabilizer Diluent (5X CS2)			
Catalog#	Size	Lot#	Expiry at 2-8°C
586	Bulk	17F16	02/2019
6169	25 mL	NA	NA
6102	100 mL	17F19	02/2019
6103	500 mL	17F18	02/2019
6104	1 L	17F17	02/2019
6105	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	62,100	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control 
Quality Assurance

Date 02/12/17
02/12/2017