

CERTIFICATE OF ANALYSIS

5X Antigen-Down HRP Conjugate Stabilizer Diluent (5X CS2)			
Catalog#	Size	Lot#	Expiry at 2-8°C
586	Bulk	14N20	06/2016
6169	25 mL	NA	NA
6102	100 mL	NA	NA
6103	500 mL	14N21 15N9	06/2016 06/2016
6104	1 L	14N22	06/2016
6105	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH at 5X	7.2-7.6	7.39	Pass
Appearance	Clear to slightly cloudy, yellow tinted liquid	Conforms	Pass
Conductivity uMHos	60,000-80,000 uMHO	76,000	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Hanson
 Tracy Hanson

Date 06/02/2015
 06/02/2015