

**CERTIFICATE OF ANALYSIS**

Page 1 of 3

**Product Description:** Kidney Injury Panel 3 (human) Kit  
**Kit Catalog Numbers:** K15189D-Series  
**Kit Lot Number:** K0040439  
**Expiration Date:** 30 Jun 2019

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		KIP-3 Control 1		KIP-3 Control 2	
		Recovery	CV	Recovery	CV	Recovery	CV
1	αGST	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	Calbindin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	Clusterin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	KIM-1	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
5	Osteoactivin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	TFF3	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
7	VEGF	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

**Acceptance Data/ Non Specific Binding Specifications:**

Spot	Name	NSB Spec
1	αGST	< 1.0%
2	Calbindin	< 1.0%
3	Clusterin	< 1.0%
4	KIM-1	< 1.0%
5	Osteoactivin	< 1.0%
6	TFF3	< 1.0%
7	VEGF	< 1.0%

## CERTIFICATE OF ANALYSIS

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	αGST	20X Calibrator	40000	pg/mL
		KIP-3 Control 1	8690	pg/mL
		KIP-3 Control 2	463	pg/mL
2	Calbindin	20X Calibrator	644000	pg/mL
		KIP-3 Control 1	93700	pg/mL
		KIP-3 Control 2	5810	pg/mL
3	Clusterin	20X Calibrator	4960000	pg/mL
		KIP-3 Control 1	696000	pg/mL
		KIP-3 Control 2	45200	pg/mL
4	KIM-1	20X Calibrator	530000	pg/mL
		KIP-3 Control 1	73200	pg/mL
		KIP-3 Control 2	4570	pg/mL
5	Osteoactivin	20X Calibrator	956000	pg/mL
		KIP-3 Control 1	142000	pg/mL
		KIP-3 Control 2	9810	pg/mL
6	TFF3	20X Calibrator	48000	pg/mL
		KIP-3 Control 1	6250	pg/mL
		KIP-3 Control 2	464	pg/mL
7	VEGF	20X Calibrator	56600	pg/mL
		KIP-3 Control 1	8690	pg/mL
		KIP-3 Control 2	599	pg/mL

Control data was generated utilizing internal reference controls.

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

**Standard Curve Data:**

αGST - Spot 1			Calbindin - Spot 2			Clusterin - Spot 3		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
2000	116,000	7.9	32200	1,180,000	1.4	248000	719,000	3.3
500	25,200	7.9	8050	247,000	2.9	62000	226,000	3.5
125	5,290	14.0	2010	37,900	4.4	15500	51,800	5.3
31.3	1,180	4.4	503	6,640	4.8	3880	13,600	5.7
7.81	393	7.2	126	1,770	2.5	969	3,600	6.3
1.95	181	4.3	31.4	655	3.7	242	1,200	2.8
0.488	133	5.6	7.86	417	5.8	60.5	510	4.4
0.00	107	8.6	0.00	325	3.9	0.00	369	5.6

**CERTIFICATE OF ANALYSIS**

Page 3 of 3

KIM-1 - Spot 4		
Conc. pg/mL	ECL Signal	%CV
26500	753,000	5.0
6630	282,000	4.6
1660	65,700	4.2
414	15,300	4.7
104	3,940	3.5
25.9	1,070	3.3
6.47	372	4.9
0.00	142	4.5

Osteoactivin - Spot 5		
Conc. pg/mL	ECL Signal	%CV
47800	224,000	4.2
12000	73,800	3.9
2990	21,300	6.1
747	5,540	3.8
187	1,840	3.3
46.7	816	2.8
11.7	614	6.2
0.00	508	7.6

TFF3 - Spot 6		
Conc. pg/mL	ECL Signal	%CV
2400	159,000	5.3
600	35,300	8.1
150	7,460	9.9
37.5	1,820	10.0
9.38	570	5.9
2.34	257	5.1
0.586	183	4.5
0.00	187	31.0

VEGF - Spot 7		
Conc. pg/mL	ECL Signal	%CV
2830	834,000	2.1
708	145,000	2.3
177	24,100	3.3
44.2	4,600	3.0
11.1	1,320	3.1
2.76	489	2.6
0.691	333	9.1
0.00	262	6.3

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Kidney Injury Panel 3 (human) Plate	Z0055975	2-8°C	30 Nov 2019
Anti-hu αGST Antibody	D0040705	2-8°C	31 Jan 2020
Anti-hu Calbindin Antibody	D0040706	2-8°C	30 Nov 2019
Anti-hu Clusterin Antibody	D0040707	2-8°C	30 Nov 2019
Anti-hu KIM-1 Antibody	D0040708	2-8°C	30 Nov 2019
Anti-hu Osteoactivin Antibody	D0040709	2-8°C	30 Nov 2019
Anti-hu TFF3 Antibody	D0040710	2-8°C	30 Nov 2019
Anti-hu VEGF Antibody	D0040711	2-8°C	31 Dec 2019
Kidney Injury Panel 3 (human) Calibrator Blend	A0040212	≤ -70°C	31 Jan 2021
Blocker A	Not Kit Specific	Not Kit Specific	N/A
Diluent 37	M0370019	≤ -10°C	30 Jun 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

**Statement:**

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	18 SEP 2018