

CERTIFICATE OF ANALYSIS

COA-06420

Product Description: Inflammation Panel 1 (rat) Kit
Kit Catalog Numbers: K15179C-1, K15179C-2, K15179C-4
Kit Lot Number: K0040216
Expiration Date: 28 Feb 2017

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Control 1		Control 2		Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Lipocalin-2	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	TSP-1	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
3	TIMP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	MCP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Lipocalin-2	0.033	80% - 120%	< 20%	27.6	75% - 125%	< 25%
2	TSP-1	29.7	80% - 120%	< 20%	766	75% - 125%	< 25%
3	TIMP-1	0.057	80% - 120%	< 20%	6.71	75% - 125%	< 25%
4	MCP-1	0.045	80% - 120%	< 20%	11.0	75% - 125%	< 25%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	Lipocalin-2	< 1.0%
2	TSP-1	< 3.0%
3	TIMP1	< 1.0%
4	MCP-1	< 1.0%

CERTIFICATE OF ANALYSIS

COA-06420

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Lipocalin-2	20X Calibrator	1160	ng/ml
		Control 1	6.15	ng/ml
		Control 2	2.66	ng/ml
		Control 3	0.0781	ng/ml
2	TSP-1	20X Calibrator	50400	ng/ml
		Control 1	266	ng/ml
		Control 2	225	ng/ml
		Control 3	59.6	ng/ml
3	TIMP-1	20X Calibrator	192	ng/ml
		Control 1	2.23	ng/ml
		Control 2	0.458	ng/ml
		Control 3	0.102	ng/ml
4	MCP-1	20X Calibrator	314	ng/ml
		Control 1	6.19	ng/ml
		Control 2	0.820	ng/ml
		Control 3	0.135	ng/ml

Controls tested may not be the same lot as shipped.

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:

Lipocalin-2 - Spot 1		
Conc. ng/ml	ECL Signal	%CV
58.0	348,000	2.0
19.3	198,000	3.8
6.44	75,000	6.1
2.15	26,700	3.1
0.716	9,070	4.9
0.239	3,160	4.5
0.0796	1,070	6.5
0.00	54	11.0

TSP-1 - Spot 2		
Conc. ng/ml	ECL Signal	%CV
2520	4,220	5.5
840	3,030	5.5
280	1,590	8.3
93.3	702	7.6
31.1	298	5.6
10.4	177	8.8
3.46	129	8.5
0.00	126	18.0

TIMP-1 - Spot 3		
Conc. ng/ml	ECL Signal	%CV
9.60	434,000	5.6
3.20	98,600	8.2
1.07	19,200	15.0
0.356	4,490	7.6
0.119	1,330	9.9
0.0395	461	4.4
0.0132	188	6.8
0.00	59	7.5

CERTIFICATE OF ANALYSIS


MCP-1 - Spot 4		
Conc. ng/ml	ECL Signal	%CV
15.7	670,000	2.0
5.23	257,000	5.2
1.74	75,500	4.3
0.581	23,700	5.0
0.194	8,410	7.6
0.0646	3,010	4.3
0.0215	1,200	5.3
0.00	316	5.0

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Inflammation Panel 1 (rat) Plate	Z0054740	2-8°C	29 Feb 2016
Anti-rat Lipocalin-2 Antibody	D0040401	2-8°C	31 Jan 2017
Anti-rat TSP-1 Antibody	D0040402	2-8°C	31 Jan 2017
Anti-rat MCP-1 Antibody	D0040403	2-8°C	31 Mar 2016
Anti-rat TIMP-1 Antibody	D0040404	2-8°C	31 Mar 2016
Inflammation Panel 1 (rat) Calibrator Blend	A0040080	≤ -70°C	31 Aug 2016
Diluent 7	M0070042	≤ -10°C	31 May 2016
Diluent 100	Y0270195	2-8°C	30 Apr 2016
Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A

Statement:

The above kit was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance		02FEB2016

CERTIFICATE OF ANALYSIS


COA-06420

Description: Inflammation Panel 1 (rat) Kit

Catalog Number(s): K15179C-1/-2/-4

Lot #: K0040216

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of Lot # K0040216 has been extended to February 28, 2017.

Function	Name	Signature	Date
Quality Assurance	Michael Grossman		01 Feb 2016

This memo supersedes all other statements of date of expiration for this kit, Lot # K0040216.