

## CERTIFICATE OF ANALYSIS

Page 1 of 3

**Product Description:** Inflammation Panel 1 (rat) Kit  
**Kit Catalog Numbers:** K15179C-Series  
**Kit Lot Number:** K0040429  
**Expiration Date:** 31 Jul 2019

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		Inflammation Panel 1 (rat) Control 1		Inflammation Panel 1 (rat) Control 2		Inflammation Panel 1 (rat) 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	0.0330	80% - 120%	< 20%	27.6	75% - 125%	< 25%
2	TSP-1	29.7	80% - 120%	< 20%	766	75% - 125%	< 25%
3	TIMP-1	0.0567	80% - 120%*	< 20%	6.71	75% - 125%	< 25%
4	MCP-1	0.0447	80% - 120%	< 20%	11.0	75% - 125%	< 25%

\* See Additional Comments.

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

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

## CERTIFICATE OF ANALYSIS

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Lipocalin-2	20X Calibrator	1032	ng/mL
		Inflammation Panel 1 (rat) Control 1	225	ng/mL
		Inflammation Panel 1 (rat) Control 2	20.4	ng/mL
		Inflammation Panel 1 (rat) Control 3	2.12	ng/mL
2	TSP-1	20X Calibrator	85900	ng/mL
		Inflammation Panel 1 (rat) Control 1	9150	ng/mL
		Inflammation Panel 1 (rat) Control 2	5500	ng/mL
		Inflammation Panel 1 (rat) Control 3	3290	ng/mL
3	TIMP-1	20X Calibrator	193	ng/mL
		Inflammation Panel 1 (rat) Control 1	66.9	ng/mL
		Inflammation Panel 1 (rat) Control 2	17.6	ng/mL
		Inflammation Panel 1 (rat) Control 3	3.65	ng/mL
4	MCP-1	20X Calibrator	292	ng/mL
		Inflammation Panel 1 (rat) Control 1	88.1	ng/mL
		Inflammation Panel 1 (rat) Control 2	18.5	ng/mL
		Inflammation Panel 1 (rat) Control 3	3.63	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			TSP-1 - Spot 2			TIMP-1 - Spot 3		
Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV
51.6	98,700	4.0	4300	17,800	4.8	9.65	239,000	4.6
17.2	56,400	5.7	1430	10,700	4.1	3.22	49,800	5.2
5.73	23,100	5.0	477	4,520	3.4	1.07	8,280	4.7
1.91	7,960	4.2	159	1,590	4.4	0.357	1,600	5.0
0.637	2,520	3.8	53.0	602	5.3	0.119	432	5.3
0.212	860	6.3	17.7	300	6.5	0.0397	160	7.4
0.0708	321	4.6	5.89	200	9.5	0.0132	85	7.3
0.00	52	9.4	0.00	163	19.0	0.00	48	9.7

**CERTIFICATE OF ANALYSIS**

MCP-1 - Spot 4		
Conc. ng/mL	ECL Signal	%CV
14.6	462,000	2.4
4.87	239,000	2.4
1.62	76,900	3.2
0.541	23,900	4.0
0.180	7,960	3.5
0.0601	2,790	4.4
0.0200	1,160	4.6
0.00	326	2.5

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Inflammation Panel 1 (rat) Plate	Z0056038	2-8°C	31 Jan 2020
Anti-rat Lipocalin-2 Antibody	D0040735	2-8°C	29 Feb 2020
Anti-rat TSP-1 Antibody	D0040770	2-8°C	30 Nov 2021
Anti-rat TIMP-1 Antibody	D0040771	2-8°C	30 Nov 2021
Anti-rat MCP-1 Antibody	D0040737	2-8°C	31 Mar 2020
Inflammation Panel 1 (rat) Calibrator Blend	A0040189	≤ -70°C	31 Dec 2019
Diluent 7	M0070052	≤ -10°C	31 Jul 2019
Diluent 100	Y0270253	2-8°C	30 Apr 2020
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

**Additional Comments:**

	Assay	Spot	Specification	Comment
1	Lipocalin-2	1	50-200%	Measured Avg Signal Ratio = 29.9%
2	TIMP-1	3	50-200%	Measured Avg Signal Ratio = 27.8%
3	TIMP-1	3	80-120%	Measured Recovery of LLOQ Conc.= 124.7%.

**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>	21 AUG 2018