

CERTIFICATE OF ANALYSIS

Page 1 of 3

Product Description: Inflammation Panel 1 (rat) Kit
Kit Catalog Numbers: K15179C-Series
Kit Lot Number: K0040473
Expiration Date: 31 Mar 2021

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%

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	1072	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	43600	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	186	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	266	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

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:

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
53.6	610,000	7.6	2180	25,700	5.7	9.32	247,000	4.8
17.9	379,000	2.3	727	11,500	7.3	3.11	53,600	5.5
5.96	165,000	3.9	242	4,170	7.6	1.04	9,470	6.8
1.99	59,100	2.5	80.7	1,430	6.3	0.345	1,860	3.3
0.662	20,700	3.9	26.9	539	7.8	0.115	564	6.4
0.221	7,110	3.1	8.97	233	4.7	0.0384	219	7.0
0.0735	2,500	3.8	2.99	121	4.9	0.0128	148	16.0
0.00	132	13.0	0.00	68	11.0	0.00	80	22.0

CERTIFICATE OF ANALYSIS


MCP-1 - Spot 4		
Conc. ng/mL	ECL Signal	%CV
13.3	396,000	2.8
4.43	166,000	4.9
1.48	49,600	7.5
0.493	15,400	3.8
0.164	5,170	4.3
0.0547	1,930	3.2
0.0182	892	2.2
0.00	358	2.2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Inflammation Panel 1 (rat) Plate	Z0056531	2-8°C	30 Nov 2021
Anti-rat Lipocalin-2 Antibody	D0040858	2-8°C	30 Nov 2023
Anti-rat TSP-1 Antibody	D0040859	2-8°C	30 Jun 2023
Anti-rat TIMP-1 Antibody	D0040861	2-8°C	30 Nov 2022
Anti-rat MCP-1 Antibody	D0040860	2-8°C	30 Nov 2022
Inflammation Panel 1 (rat) Calibrator Blend	A0040220	≤ -70°C	31 Mar 2021
Diluent 100	Not Kit Specific	Not Kit Specific	N/A
Diluent 7	M0070060	≤ -10°C	31 Oct 2021
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
Review/Approval	Karen Hamilla	Quality		24 JAN 2020