

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Inflammation Panel 1 (rat) Kit
Kit Catalog Numbers: K15179C-Series
Kit Lot Number: K0040445
Expiration Date: 31 Jan 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%

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

Controls data was generated using internal reference controls.

CERTIFICATE OF ANALYSIS

Page 2 of 2

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	500,000	6.7	2180	43,600	5.4	9.32	256,000	6.0
17.9	329,000	7.0	727	18,600	7.0	3.11	53,300	10.0
5.96	142,000	6.5	242	6,180	12.0	1.04	8,370	11.0
1.99	50,500	6.2	80.7	1,940	13.0	0.345	1,540	11.0
0.662	16,600	7.2	26.9	703	7.2	0.115	466	8.6
0.221	5,990	5.9	8.97	318	8.1	0.0384	185	11.0
0.0735	2,090	5.3	2.99	170	5.2	0.0128	108	5.2
0.00	75	12.0	0.00	106	6.5	0.00	85	20.0


MCP-1 - Spot 4		
Conc. ng/mL	ECL Signal	%CV
13.3	241,000	2.6
4.43	124,000	4.1
1.48	40,100	4.5
0.493	12,300	6.9
0.164	4,340	2.7
0.0547	1,580	5.3
0.0182	740	3.4
0.00	261	3.1

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Inflammation Panel 1 (rat) Plate	Z0056360	2-8°C	28 Feb 2021
Anti-rat Lipocalin-2 Antibody	D0040803	2-8°C	30 Sep 2022
Anti-rat TSP-1 Antibody	D0040804	2-8°C	31 Aug 2022
Anti-rat TIMP-1 Antibody	D0040806	2-8°C	30 Jun 2022
Anti-rat MCP-1 Antibody	D0040805	2-8°C	28 Feb 2022
Inflammation Panel 1 (rat) Calibrator Blend	A0040220	≤ -70°C	31 Mar 2021
Diluent 7	M0070056	≤ -10°C	31 Jan 2021
Diluent 100	Y0270260	2-8°C	30 Jun 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	Dawn Stollar	Quality		29 MAY 2019