

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Inflammation Panel 3 (cyno) Kit
Kit Catalog Numbers: K15191D-Series
Kit Lot Number: K0040490
Expiration Date: 30 Jun 2022

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Inflammation Panel 3 Control 1		Inflammation Panel 3 Control 2		Inflammation Panel 3 Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	MCP-1	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
2	NGAL	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
3	TIMP-1	80% - 120%	< 20%	75% - 125%	< 25%	80% - 125%	< 25%	N/A	N/A

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	MCP-1	< 1.0%
2	NGAL	< 1.0%
3	TIMP-1	< 1.0%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	MCP-1	20X Calibrator	18	ng/mL
		Inflammation Panel 3 Control 1	4.29	ng/mL
		Inflammation Panel 3 Control 2	2.13	ng/mL
		Inflammation Panel 3 Control 3	1.12	ng/mL
2	NGAL	20X Calibrator	504	ng/mL
		Inflammation Panel 3 Control 1	87.4	ng/mL
		Inflammation Panel 3 Control 2	39.7	ng/mL
		Inflammation Panel 3 Control 3	21.6	ng/mL
3	TIMP-1	20X Calibrator	1076	ng/mL
		Inflammation Panel 3 Control 1	286	ng/mL
		Inflammation Panel 3 Control 2	91.1	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

CERTIFICATE OF ANALYSIS

Standard Curve Data:


MCP-1 - Spot 1			NGAL - Spot 2			TIMP-1 - Spot 3		
Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV
0.901	552,000	1.6	25.2	208,000	1.9	53.8	236,000	4.2
0.300	370,000	1.7	8.40	95,100	1.4	17.9	73,700	2.8
0.100	152,000	1.6	2.80	34,800	1.7	5.98	10,300	3.8
0.0334	51,600	2.5	0.933	12,500	2.0	1.99	1,370	3.0
0.0111	14,800	1.9	0.311	4,290	2.9	0.664	333	2.8
0.00371	4,110	1.6	0.104	1,590	1.9	0.221	129	3.5
0.00124	1,150	1.7	0.0346	624	3.0	0.0738	76	5.7
0.00	51	12.0	0.00	109	14.0	0.00	50	9.6

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Inflammation Panel 3 (Cyno) Plate	Z0056668	2-8°C	31 Jul 2022
Anti-hu MCP-1 Antibody	D0040904	2-8°C	31 Aug 2023
Anti-hu NGAL Antibody	D0040906	2-8°C	31 Aug 2023
Anti-hu TIMP-1 Antibody	D0040905	2-8°C	30 Sep 2023
Inflammation Panel 3(Cyno) Calibrator Blend	A0040248	≤ -70°C	31 Oct 2022
Blocker A	Not Kit Specific	Not Kit Specific	N/A
Diluent 8	M0080036	≤ -10°C	30 Sep 2022
Diluent 11	M0110044	≤ -10°C	30 Jun 2022
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		28 AUG 2020