

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Inflammation Panel 3 (cyno) Kit

Kit Catalog Numbers: K15191D-Series

Kit Lot Number: K0040428

Expiration Date: 31 Oct 2020

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Control 1		Control 2		Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	MCP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	NGAL	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	TIMP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

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	17.4	ng/mL
		Control 1	0.291	ng/mL
		Control 2	0.0969	ng/mL
		Control 3	0.0323	ng/mL
2	NGAL	20X Calibrator	522	ng/mL
		Control 1	8.7	ng/mL
		Control 2	2.9	ng/mL
		Control 3	0.967	ng/mL
3	TIMP-1	20X Calibrator	1038	ng/mL
		Control 1	17.3	ng/mL
		Control 2	5.77	ng/mL
		Control 3	1.92	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
0.872	733,000	1.7	26.1	239,000	2.3	51.9	288,000	2.3
0.291	400,000	2.3	8.70	109,000	2.0	17.3	102,000	1.8
0.0969	153,000	1.7	2.90	40,200	3.1	5.77	16,400	5.5
0.0323	50,600	1.0	0.967	13,900	3.9	1.92	2,200	3.7
0.0108	14,400	1.6	0.322	4,710	3.7	0.641	500	3.7
0.00359	4,200	2.9	0.107	1,720	3.9	0.214	208	2.5
0.00120	1,310	2.4	0.0358	676	1.4	0.0712	134	4.4
0.00	96	6.1	0.00	166	13.0	0.00	99	6.1

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Inflammation Panel 3 (cyno) Plate	Z0056267	2-8°C	31 Oct 2020
Anti-hu MCP-1 Antibody	D0040767	2-8°C	31 Oct 2021
Anti-hu NGAL Antibody	D0040769	2-8°C	31 Jan 2022
Anti-hu TIMP-1 Antibody	D0040768	2-8°C	31 Oct 2021
Inflammation Panel 3 (cyno) Calibrator Blend	A0040214	≤ -70°C	30 Nov 2020
Blocker A	Not Kit Specific	Not Kit Specific	N/A
Diluent 8	M0080033	≤ -10°C	30 Nov 2020
Diluent 11	M0110035	≤ -10°C	31 Dec 2020
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 Assurance		30 OCT 2018