

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Inflammation Panel 3 (cyno) Kit  
**Kit Lot Number:** K0040507  
**Expiration Date:** 31 Jan 2022  
**Kit Catalog Numbers:** K15191D-Series

**Precision and Accuracy 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%	75% - 125%	< 25%	N/A	N/A

**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 Calibrator:**


Spot	Name	Conc.	Units
1	MCP-1	18	ng/mL
2	NGAL	504	ng/mL
3	TIMP-1	1076	ng/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Inflammation Panel 3 (cyno) Plate	Z0056784	2-8°C	28 Feb 2023
Anti-hu MCP-1 Antibody	D0040935	2-8°C	31 Mar 2024
Anti-hu NGAL Antibody	D0040937	2-8°C	29 Feb 2024
Anti-hu TIMP-1 Antibody	D0040936	2-8°C	30 Apr 2024
Inflammation Panel 3 (cyno) Calibrator Blend	A0040248	≤ -70°C	31 Oct 2022
Diluent 8	M0080038	≤ -10°C	31 Dec 2022
Diluent 11	M0110041	≤ -10°C	31 Jan 2022

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		21 DEC 2021