

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** Human TIMP-1 Kit  
**Kit Catalog Numbers:** K151JFC-Series  
**Kit Lot Number:** K0040464  
**Expiration Date:** 30 Jun 2021

**Acceptance Data/ Percent Recovery Specifications:**

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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TIMP-1	494	80% - 120%	< 20%	32000	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	TIMP-1	10X Calibrator	406000	pg/mL
		Human TIMP-1 Control 1	835000	pg/mL
		Human TIMP-1 Control 2	249000	pg/mL
		Human TIMP-1 Control 3	83000	pg/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:**

TIMP-1 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
40600	1,260,000	1.7
13500	397,000	3.9
4510	82,300	2.9
1500	13,100	4.4
501	2,280	5.1
167	558	6.7
55.7	236	28.0
0.00	122	70.0


**CERTIFICATE OF ANALYSIS**

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human TIMP-1 Plate	Z0056492	2-8°C	30 Sep 2021
Anti-Hu TIMP-1 Antibody	D0040835	2-8°C	30 Nov 2022
Human TIMP-1 Calibrator	A0040224	≤ -70°C	30 Jun 2021
Diluent 2	M0020106	≤ -10°C	30 Sep 2021
Diluent 3	M0030089	≤ -10°C	30 Sep 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
<b>Review/Approval</b>	Karen Hamilla	Quality		24 SEP 2019