

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Human TIMP-1 Kit
Kit Catalog Numbers: K151JFC-Series
Kit Lot Number: K0040482
Expiration Date: 31 Jan 2022

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	419000	pg/mL
		Human TIMP-1 Control 1	93500	pg/mL
		Human TIMP-1 Control 2	43500	pg/mL
		Human TIMP-1 Control 3	17900	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
41900	1,020,000	2.5
14000	304,000	3.7
4660	63,600	6.1
1550	10,400	6.4
517	1,780	6.5
172	410	6.5
57.5	148	1.9
0.00	62	5.9

Components:


CERTIFICATE OF ANALYSIS

Page 2 of 2

Description	Lot Number	Storage Temperature	Expiration Date
Human TIMP-1 Plate	Z0056607	2-8°C	31 Mar 2022
Anti-hu TIMP-1 Antibody	D0040881	2-8°C	30 Sep 2023
Human TIMP-1 Calibrator	A0040241	≤ -70°C	30 Apr 2022
Diluent 2	M0020107	≤ -10°C	31 Jan 2022
Diluent 3	M0030097	≤ -10°C	31 Aug 2022

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		03 FEB 2021