

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Human TIMP-1 Kit
Kit Catalog Numbers: K151JFC-Series
Kit Lot Number: K0040451
Expiration Date: 30 Apr 2020

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
		Control 1	835000	pg/mL
		Control 2	249000	pg/mL
		Control 3	83000	pg/mL

Controls 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	1,290,000	1.7
13500	407,000	2.1
4510	80,900	4.9
1500	12,700	9.6
501	2,010	4.9
167	435	5.4
55.7	169	3.7
0.00	86	4.9

CERTIFICATE OF ANALYSIS


Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TIMP-1 Plate	Z0056398	2-8°C	31 Mar 2021
Human TIMP-1 Antibody	D0040812	2-8°C	30 Apr 2022
Human TIMP-1 Calibrator	A0040224	≤ -70°C	30 Jun 2021
Diluent 2	M0020102	≤ -10°C	28 Feb 2021
Diluent 3	M0030077	≤ -10°C	30 Apr 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		18 MAR 2019