

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Human TIMP-1 Kit

Kit Catalog Numbers: K151JFC-Series

Kit Lot Number: K0040413

Expiration Date: 30 Sep 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Control 1		Control 2		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	410000	pg/mL
		Control 1	835000	pg/mL
		Control 2	249000	pg/mL
		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
41000	1,420,000	1.1
13700	488,000	5.4
4560	98,000	4.0
1520	15,500	4.9
506	2,630	5.8
169	585	4.6
56.2	200	4.6
0.00	82	5.8

CERTIFICATE OF ANALYSIS


Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TIMP-1 Plate	Z0056071	2-8°C	29 Feb 2020
Anti-hu TIMP-1 Antibody	D0040743	2-8°C	29 Feb 2020
Human TIMP-1 Calibrator	A0040204	≤ -70°C	29 Feb 2020
Diluent 2	M0020096	≤ -10°C	30 Nov 2019
Diluent 3	M0030076	≤ -10°C	30 Sep 2019
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		07 MAR 2019