

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Human TIMP-1 Kit
Kit Catalog Numbers: K151JFC-series
Kit Lot Number: K0040433
Expiration Date: 30 Sep 2019

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	350000	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
35000	1,160,000	3.6
11700	340,000	5.9
3890	64,000	6.9
1300	10,900	4.4
432	1,960	4.1
144	493	5.5
48.0	191	5.0
0.00	78	7.3

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TIMP-1 Plate	Z0056298	2-8°C	31 Dec 2020
Anti-hu TIMP-1 Antibody	D0040780	2-8°C	31 Dec 2021
Human TIMP-1 Calibrator	A0040217	≤ -70°C	31 Dec 2020
Diluent 2	M0020101	≤ -10°C	31 Dec 2020
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	<i>Dawn Stollar</i>	28 SEP 2018