



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-05040

Product Description: Human TIMP-1 Kit

Kit Catalog Numbers: K151JFC-1, K151JFC-2, K151JFC-4

Kit Lot Number: K0040277

Expiration Date: 31 DEC 2016

Acceptance Data:

Metric		Specifications	Assay
			Human TIMP-1
% Recovery Concentration	Calibrators	80% - 120%	PASS
	Controls	80% - 120%	PASS
	ULOQ	80% - 120%	PASS
	LLOQ	80% - 120%	PASS
% Recovery CV	Calibrators	< 20%	PASS
	Controls	< 20%	PASS
	ULOQ	< 20%	PASS
	LLOQ	< 20%	PASS

Concentration of Materials Tested:

Units = pg/mL	Assay
Human TIMP-1	
10X Calibrator ¹	382 000
Control 1	715 000
Control 2	211 500
Control 3	70 500
ULOQ	32 000
LLOQ	494

Standard Curve Data from Kit QC:

Human TIMP-1		
Conc., (pg/mL)	Average Signal	%CV
38 200	1 343 441	4.5
12 733	439 574	6.4
4244	101 527	7.5
1415	17 359	9.1
472	3521	9.6
157	707	7.0
52.4	223	8.4
0	97	25.9

¹ Concentration provided in this table refer to the stock concentration of the calibrator. For top of the standard curve concentration, please refer to the Standard Curve Data on page 1.




CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-05040

Human TIMP-1 Kit Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human TIMP-1 Plate	Z0055206	2-8°C	30 APR 2017
Anti-hu TIMP-1 Antibody	D0040496	2-8°C	30 APR 2017
Human TIMP-1 Calibrator	A0040093	≤ -70°C	31 DEC 2016
Diluent 2	M0020085	≤ -10°C	31 JAN 2017
Diluent 3	M0030063	≤ -10°C	30 APR 2017
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A

Statement: The above kit lot was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dana Short	Quality Control		19 DEC 2014