

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human TIMP-1 Kit
Kit Lot Number: K0040514
Expiration Date: 31 Jul 2023
Kit Catalog Numbers: K151JFC-series

Precision and Accuracy Specifications:

		Calibrator		TIMP-1 Control 1		TIMP-1 Control 2		TIMP-1 Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	TIMP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	TIMP-1	494	80% - 120%	< 20%	32000	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	TIMP-1	468000	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TIMP-1 Plate	Z0056869	2-8°C	31 Jul 2023
Anti-hu TIMP-1 Antibody	D0040948	2-8°C	28 Feb 2025
Human TIMP-1 Calibrator	A0040260	≤ -70°C	30 Sep 2023
Diluent 2	M0020117	≤ -10°C	31 Oct 2023
Diluent 3	M0030107	≤ -10°C	30 Sep 2023

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 OCT 2021