

**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** Human TIMP-1 Kit  
**Kit Lot Number:** K0040535  
**Expiration Date:** 31 Oct 2024  
**Kit Catalog Numbers:** K151JFC-Series

**Precision and Accuracy 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%	180% - 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 Calibrator:**


Spot	Name	Conc.	Units
1	TIMP-1	420000	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human TIMP-1 Plate	Z0057224	2-8°C	31 Jan 2025
Anti-hu TIMP-1 Antibody	D0040999	2-8°C	31 May 2026
Human TIMP-1 Calibrator	A0040265	≤ -70°C	31 Oct 2024
Diluent 2	M0020123	≤ -10°C	28 Feb 2025
Diluent 3	M0030118	≤ -10°C	31 Mar 2025

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		07 DEC 2022