



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human TIMP-1 Calibrator

Catalog Number: C01JF-2

Lot Number: A0040260

Expiration Date: 30 SEP 2023


Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
TIMP-1	468000	pg/mL

Dilution to Top of Curve: 10X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		05 OCT 2021