



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human TIMP-1 Calibrator

Catalog Number: C01JF-2

Lot Number: A0040224

Expiration Date: 30 JUN 2021

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

Concentration of Materials Tested:

	Assay
	TIMP-1
Calibrator (pg/mL)	406000

Concentration: 10X

Statement: This 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>	04 MAR 2019