



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human Tissue Culture Calibrator 2

Catalog Number: C0105-2

Lot Number: A0032444

Expiration Date: 31 Mar 2022

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

Concentration Assignments:

	Nominal Concentration		Measured Concentration	
	Stock Concentration ($\mu\text{g/mL}$)	Top of Curve Concentration (pg/mL)	Stock Concentration ($\mu\text{g/mL}$)	Top of Curve Concentration (pg/mL)
IL-1 β	1.00	10000	1.02	10200
IL-2	1.00	10000	1.14	11400
IL-4	1.00	10000	1.05	10500
IL-5	1.00	10000	1.01	10100
IL-6	1.00	10000	0.98	9750
IL-8	1.00	10000	0.91	9090
IL-10	1.00	10000	0.91	9070
IL-12p70	1.00	10000	1.48	14750
IL-13	1.00	10000	0.97	9730
GM-CSF	1.00	10000	0.95	9540
TNF- α	1.00	10000	1.03	10280

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

	Name	Function	Signature	Date
Review/Approval	Xiomara Talavera	Quality		03 JAN 2020