



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human Tissue Culture Calibrator 2

Catalog Number: C0105-2

Lot Number: A0032399

Expiration Date: 28 FEB 2021

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	10,000	0.92	9,230
IL-2	1.00	10,000	0.95	9,490
IL-4	1.00	10,000	0.90	8,960
IL-5	1.00	10,000	0.95	9,490
IL-6	1.00	10,000	0.86	8,630
IL-8	1.00	10,000	0.88	8,790
IL-10	1.00	10,000	0.94	9,380
IL-12p70	1.00	10,000	0.92	9,200
IL-13	1.00	10,000	1.05	10,500
GM-CSF	1.00	10,000	0.96	9,570
TNF- α	1.00	10,000	1.00	9,980

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 JAN 2019