



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human Tissue Culture Calibrator 2

Catalog Number: C0105-2

Lot Number: A0032297

Expiration Date: 31 Oct 2019

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	0.99	9 930
IL-2	1.00	10,000	1.03	10 300
IL-4	1.00	10,000	1.02	10 200
IL-5	1.00	10,000	1.03	10 300
IL-6	N/A	N/A	N/A	N/A
IL-8	1.00	10,000	0.99	9 930
IL-10	1.00	10,000	1.02	10 200
IL-12p70	1.00	10,000	1.00	10 040
IL-13	1.00	10,000	0.96	9 600
GM-CSF	1.00	10,000	1.04	10 400
TNF- α	1.00	10,000	0.97	9 710

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>	05 JUL 2017