



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human Tissue Culture Calibrator 2

Catalog Number: C0105-2

Lot Number: A0032316

Expiration Date: 30 Nov 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	0.91	9 100
IL-2	1.00	10 000	1.08	10 800
IL-4	1.00	10 000	1.01	10 100
IL-5	1.00	10 000	1.03	10 300
IL-6	1.00	10 000	0.97	9 670
IL-8	1.00	10 000	0.99	9 880
IL-10	1.00	10 000	1.03	10 300
IL-12p70	1.00	10 000	1.02	10 200
IL-13	1.00	10 000	0.94	9 400
GM-CSF	1.00	10 000	1.05	10 500
TNF- α	1.00	10 000	1.01	10 080

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

	Name	Function	Signature	Date
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	05JUL2017