



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human Tissue Culture Calibrator 2

Catalog Number: 3-3004-020304-A

Lot Number: A0032285

Expiration Date: 30 Apr 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	1.06	10,600
IL-2	1.00	10,000	1.16	11,600
IL-4	1.00	10,000	1.05	10,500
IL-5	1.00	10,000	1.08	10,800
IL-6	1.00	10,000	1.08	10,800
IL-8	1.00	10,000	1.00	10,030
IL-10	1.00	10,000	1.03	10,300
IL-12p70	1.00	10,000	1.06	10,600
IL-13	1.00	10,000	0.98	9,800
GM-CSF	1.00	10,000	1.00	10,000
TNF- α	1.00	10,000	1.02	10,200

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>	08 DEC 2016