



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA - 07285

Product Description: Human Tissue Culture Calibrator 2

Catalog Number: C01XX-2

Lot Number: A0032258

Expiration Date: 31 JAN 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	1.03	10,300
IL-2	1.00	10,000	1.12	11,200
IL-4	1.00	10,000	1.00	9,990
IL-5	1.00	10,000	1.06	10,600
IL-6	1.00	10,000	1.01	10,100
IL-8	1.00	10,000	0.96	9,610
IL-10	1.00	10,000	0.99	9,930
IL-12p70	1.00	10,000	0.99	9,850
IL-13	1.00	10,000	0.91	9,070
GM-CSF	1.00	10,000	0.97	9,730
TNF- α	1.00	10,000	0.97	9,720

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. 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>	12 SEP 2016