



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Human Tissue Culture Calibrator 2

**Catalog Number:** C0105-2

**Lot Number:** A0032347

**Expiration Date:** 31 Jan 2020

**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

#### Concentration Assignments:

	Nominal Concentration		Measured Concentration	
	Stock Concentration ( $\mu\text{g/mL}$ )	Top of Curve Concentration ( $\text{pg/mL}$ )	Stock Concentration ( $\mu\text{g/mL}$ )	Top of Curve Concentration ( $\text{pg/mL}$ )
IL-1 $\beta$	1.00	10,000	1.00	10,000
IL-2	1.00	10,000	0.97	9,720
IL-4	1.00	10,000	0.98	9,780
IL-5	1.00	10,000	0.98	9,810
IL-6	1.00	10,000	0.96	9,570
IL-8	1.00	10,000	0.97	9,740
IL-10	1.00	10,000	1.04	10,350
IL-12p70	1.00	10,000	0.99	9,920
IL-13	1.00	10,000	0.96	9,630
GM-CSF	1.00	10,000	1.00	9,980
TNF- $\alpha$	1.00	10,000	0.95	9,550

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	08SEP2017