



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0105-2

**Lot Number:** A0032473

**Expiration Date:** 31 JUL 2023

**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-1b	1.00	10000	0.95	9460
IL-2	1.00	10000	1.09	10850
IL-4	1.00	10000	1.06	10600
IL-5	1.00	10000	1.01	10100
IL-6	1.00	10000	0.86	8630
IL-8	1.00	10000	0.91	9110
IL-10	1.00	10000	0.97	9660
IL-12p70	1.00	10000	1.00	10010
IL-13	1.00	10000	1.10	11000
GM-CSF	1.00	10000	0.91	9070
TNF-a	1.00	10000	0.94	9440

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		01 APR 2021