



### CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0105-2

**Lot Number:** A0032493

**Expiration Date:** 31 DEC 2024

**Storage Temperature:** ≤ -70°C

**Concentration Assignments:**

	Nominal Concentration		Measured Concentration	
	Stock Concentration (µg/mL)	Top of Curve Concentration (pg/mL)	Stock Concentration (µg/mL)	Top of Curve Concentration (pg/mL)
IL-1β	1.00	10000	0.911	9110
IL-2	1.00	10000	1.12	11200
IL-4	1.00	10000	1.11	11100
IL-5	1.00	10000	0.999	9990
IL-6	1.00	10000	0.972	9720
IL-8	1.00	10000	1.005	10050
IL-10	1.00	10000	1.045	10450
IL-12p70	1.00	10000	1.085	10850
IL-13	1.00	10000	1.05	10500
GM-CSF	1.00	10000	1.05	10500
TNF-α	1.00	10000	0.948	9480

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Xiomara Talavera	Quality		27 JUL 2022