



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human Tissue Culture Calibrator 2

Catalog Number: C0105-2

Lot Number: A0032460

Expiration Date: 31 OCT 2022

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 | 1.08 | 10,800 |
| IL-5 | 1.00 | 10,000 | 1.09 | 10,900 |
| IL-6 | 1.00 | 10,000 | 0.99 | 9,880 |
| IL-8 | 1.00 | 10,000 | 0.96 | 9,620 |
| IL-10 | 1.00 | 10,000 | 1.07 | 10,650 |
| IL-12p70 | 1.00 | 10,000 | 1.10 | 11,000 |
| IL-13 | 1.00 | 10,000 | 1.17 | 11,700 |
| GM-CSF | 1.00 | 10,000 | 1.03 | 10,300 |
| TNF- α | 1.00 | 10,000 | 1.03 | 10,300 |

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

| | Name | Function | Signature | Date |
|-----------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality | | 07 MAY 2020 |