



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human IFN- γ Calibrator

Catalog Number: C01AE-2

Lot Number: A0032403

Expiration Date: 31 JAN 2021

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

Concentration Assignments:

| | Nominal Concentration | | Measured Concentration | |
|---------------|--|--|--|--|
| | Stock Concentration ($\mu\text{g}/\text{mL}$) | Top of Curve Concentration (pg/mL) | Stock Concentration ($\mu\text{g}/\text{mL}$) | Top of Curve Concentration (pg/mL) |
| IFN- γ | 1.00 | 10,000 | 1.16 | 11,600 |

Statement: 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> | 07 SEP 2018 |