



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 20  
**Catalog Number:** C0328-2  
**Lot Number:** A00U0294  
**Expiration Date:** 31 OCT 2027  
**Storage Temperature:** 2-8°C  
**Intended Assay Species:** Human

**Concentration Assignment:**

| Analyte          | Amount of Analyte in Calibrator | Unit of Measure | Stock Concentration of Reconstituted Calibrator | Unit of Measure |
|------------------|---------------------------------|-----------------|---|-----------------|
| CD276/B7-H3      | 101250                          | pg              | 405000  | pg/mL           |
| CD20             | 196000                          | pg              | 784000  | pg/mL           |
| CD27             | 9125                            | pg              | 36500   | pg/mL           |
| CD28             | 337500                          | pg              | 1350000   | pg/mL           |
| CTLA-4           | 3475                            | pg              | 13900   | pg/mL           |
| GITRL/TNFSF18    | 2700                            | pg              | 10800   | pg/mL           |
| OX40/TNFRSF4     | 4375                            | pg              | 17500   | pg/mL           |
| PD-1 (epitope 1) | 2525                            | pg              | 10100   | pg/mL           |
| PD-1 (epitope 2) | 2525                            | pg              | 10100   | pg/mL           |
| TIGIT            | 7975                            | pg              | 31900   | pg/mL           |
| TLR1             | 91000                           | pg              | 364000  | pg/mL           |

**Reconstitution Volume:** 250 µL

To calculate the **Calibrator Standard 1 (top of curve) concentration**, please refer to the dilution-factor used in the **Prepare Calibrator Standards** section of the respective product insert. Divide the stock concentration of the reconstituted calibrator by this dilution factor to generate the working calibrator standard 1 concentration.

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. 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  | <i>X Talavera</i> | 29 NOV 2022 |