



CERTIFICATE OF ANALYSIS

Product Description: Human Proinsulin Calibrator
Catalog Number: C016M-2
Lot Number: A00U0304
Expiration Date: 30 NOV 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 |
|------------|---------------------------------|-----------------|---|-----------------|
| Proinsulin | 0.2365 | pmol | 946 | pM |

Reconstitution Volume: 250 μ L

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 |
|------------------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality | | 04 JAN 2023 |