



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 15
Catalog Number: C0273-2
Lot Number: A00U0430
Expiration Date: 30 APR 2030
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
Insulin	1900	μIU	7600	μIU/mL
C-peptide	129000	pg	516000	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 is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		05 JUN 2025