



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 15

**Catalog Number:** C0273-2

**Lot Number:** A00U0068

**Expiration Date:** 31 JUL 2023

**Storage Condition:** 2-8°C

Analyte Description	Amount of Analyte in Calibrator	Units	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL)	Units
hu Insulin	1 775	µIU	7 100	µIU/mL
hu C-Peptide	71 500	pg	286 000	pg/mL

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.

**Statement:** The above calibrator lot is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	30 AUG 2018