



EXTENSION MEMO


Page 1 of 1

Product Description(s): Human GIP (active) Calibrator

Catalog Number(s) / Size(s): C016N-2

Lot Number: A00U0112

Based on additional testing, the original date of expiration for this calibrator has been extended by 12 months. The expiration date of A00U0112 has been extended to 31 MAY 2025.

Function	Name	Signature	Date
Quality	Karen Hamilla		11 APR 2024

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human GIP (active) Calibrator

Catalog Number: C016N-2

Lot Number: A00U0112

Expiration Date: 31 MAY 2024

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
GIP (active)	6 975	pg	27 900	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		30 SEP 2019