



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 2

Catalog Number: C0061-2

Lot Number: A00U0049

Expiration Date: 31 DEC 2022

Storage Condition: 2-8°C

Analyte Description	Amount of Analyte in Calibrator (pg)	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL) (pg/mL)	Concentration of Standard 1 (after 5-fold Dilution) (pg/mL)
hu Eotaxin	8 000	32 000	6 400
hu Eotaxin-3	25 750	103 000	20 600
hu IP-10	12 875	51 500	10 300
hu MCP-1	7 763	31 050	6 210
hu MCP-4	5 338	21 350	4 270
hu MDC	26 250	105 000	21 000
hu MIP-1 α	7 750	31 000	6 200
hu MIP-1 β	2 513	10 050	2 010
hu TARC	3 925	15 700	3 140

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	Dawn Stollar