



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 2

**Catalog Number:** C0061-2

**Lot Number:** A00U0011

**Expiration Date:** 31 JUL 2020

**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)
Human Eotaxin	7 488	29 950	5 990
Human Eotaxin-3	24 625	98 500	19 700
Human IP-10	11 300	45 200	9 040
Human MCP-1	8 188	32 750	6 550
Human MCP-4	5 663	22 650	4 530
Human MDC	25 125	100 500	20 100
Human MIP-1 $\alpha$	7 675	30 700	6 140
Human MIP-1 $\beta$	2 675	10 700	2 140
Human TARC	3 713	14 850	2 970

**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>	17 JUL 2017