



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05542

Product Description: Calibrator 2
Catalog Number: C0061-2
Lot Number: A00U0003
Expiration Date: 31 MAR 2020
Storage Condition: 2-8°C

Mass of Materials Tested:

	Assay								
	hu Eotaxin	hu Eotaxin-3	hu IP-10	hu MCP-1	hu MCP-4	hu MDC	hu MIP-1 α	hu MIP-1 β	hu TARC
Calibrator (pg)	6050	26 700	7500	8200	4755	25 150	5250	2045	2795

Calculated Concentration of Reconstituted Materials:

Reconstitution volume = 250 μ L

	Assay								
	hu Eotaxin	hu Eotaxin-3	hu IP-10	hu MCP-1	hu MCP-4	hu MDC	hu MIP-1 α	hu MIP-1 β	hu TARC
Calibrator Conc. (pg/mL)	24 200	106 800	30 000	32 800	19 020	100 600	21 000	8180	11 180

Statement: The above calibrator lot was tested using MSD procedures and conforms to MSD's specifications. It is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Anastasia Belilovets	Quality Assurance		24 Jul 2015