



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05069

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0001
Expiration Date: 31 OCT 2019
Storage Condition: 2-8°C

Mass of Materials Tested:

	Assay									
	hu GM-CSF	hu IFN- γ	hu IL-1 β	hu IL-2	hu IL-4	hu IL-5	hu IL-6	hu IL-8	hu IL-10	hu IL-12p70
Calibrator (pg)	11 800	21 250	4775	2420	2580	5050	2475	2760	4645	6650

	Assay			
	hu IL-13	hu IL-17A	hu TNF- α	hu VEGF-A
Calibrator (pg)	2400	29 200	4565	6150

Calculated Concentration of Reconstituted Materials:

Reconstitution volume = 250 μ L

	Assay									
	hu GM-CSF	hu IFN- γ	hu IL-1 β	hu IL-2	hu IL-4	hu IL-5	hu IL-6	hu IL-8	hu IL-10	hu IL-12p70
Calibrator Conc. (pg/mL)	47 200	85 000	19 100	9680	10 320	20 200	9900	11 040	18 580	26 600

	Assay			
	hu IL-13	hu IL-17A	hu TNF- α	hu VEGF-A
Calibrator Conc. (pg/mL)	9600	116 800	18 260	24 600

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		10 Jul 2015