



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05561

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0010
Expiration Date: 30 JUN 2020
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)	12 250	21 775	4875	2550	2483	5200	2650	2875	4875	7000

	Assay			
	hu IL-13	hu IL-17A	hu TNF- α	hu VEGF-A
Calibrator (pg)	2458	30 000	4275	5825

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)	49 000	87 100	19 500	10 200	9932	20 800	10 600	11 500	19 500	28 000

	Assay			
	hu IL-13	hu IL-17A	hu TNF- α	hu VEGF-A
Calibrator Conc. (pg/mL)	9832	120 000	17 100	23 300

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		09 Jul 2015