



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07195

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0021
Expiration Date: 31 MAY 2021
Storage Condition: 2-8°C

Mass of Materials Tested:

	Assay						
	hu IFN- γ	hu IL-1 β	hu IL-2	hu IL-4	hu IL-6	hu IL-8	hu IL-10
Calibrator (pg)	21,225	4,825	2,453	2,600	2,625	2,775	4,900

	Assay						
	hu IL-12p70	hu IL-13	hu TNF- α	hu IL-5	hu GM-CSF	hu IL-17A	hu VEGF-A
Calibrator (pg)	6,975	2,413	4,450	4,975	12,125	30,250	5,825

Calculated Concentration of Reconstituted Materials:

Reconstitution volume = 250 μ L

	Assay						
	hu IFN- γ	hu IL-1 β	hu IL-2	hu IL-4	hu IL-6	hu IL-8	hu IL-10
Calibrator (pg)	84,900	19,300	9,810	10,400	10,500	11,100	19,600

	Assay						
	hu IL-12p70	hu IL-13	hu TNF- α	hu IL-5	hu GM-CSF	hu IL-17A	hu VEGF-A
Calibrator (pg)	27,900	9,650	17,800	19,900	48,500	121,000	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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	12 AUG 2016