



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07303

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0026
Expiration Date: 30 SEP 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)	23,375	5,000	2,500	2,725	2,675	2,875	4,925

	Assay						
	hu IL-12p70	hu IL-13	hu TNF- α	hu IL-5	hu GM-CSF	hu IL-17A	hu VEGF-A
Calibrator (pg)	7,025	2,405	4,425	5,175	12,375	30,750	6,475

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/mL)	93,500	20,000	10,000	10,900	10,700	11,500	19,700

	Assay						
	hu IL-12p70	hu IL-13	hu TNF- α	hu IL-5	hu GM-CSF	hu IL-17A	hu VEGF-A
Calibrator (pg/mL)	28,100	9,620	17,700	20,700	49,500	123,000	25,900

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	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	20 SEP 2016