



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07231

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0022
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)	27,000	5,150	2,550	2,625	2,725	2,875	5,175

	Assay						
	hu IL-12p70	hu IL-13	hu TNF- α	hu IL-5	hu GM-CSF	hu IL-17A	hu VEGF-A
Calibrator (pg)	7,350	2,458	4,575	5,225	12,675	30,250	7,275

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)	108,000	20,600	10,200	10,500	10,900	11,500	20,700

	Assay						
	hu IL-12p70	hu IL-13	hu TNF- α	hu IL-5	hu GM-CSF	hu IL-17A	hu VEGF-A
Calibrator (pg)	29,400	9,830	18,300	20,900	50,700	121,000	29,100

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>	19 AUG 2016