



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05543

Product Description: Calibrator 3
Catalog Number: C0062-2
Lot Number: A00U0004
Expiration Date: 31 MAR 2020
Storage Condition: 2-8°C

Mass of Materials Tested:

	Assay									
	hu IL-1α	hu IL-7	hu IL-12/23p40	hu IL-15	hu IL-16	hu TNF-β	hu G-CSF	hu IFN-α2a	hu IL-18	hu TPO
Calibrator (pg)	6350	8800	26 250	3785	26 850	5400	25 450	53 000	52 500	50 500

Calculated Concentration of Reconstituted Materials:

Reconstitution volume = 250 µL

	Assay									
	hu IL-1α	hu IL-7	hu IL-12/23p40	hu IL-15	hu IL-16	hu TNF-β	hu G-CSF	hu IFN-α2a	hu IL-18	hu TPO
Calibrator Conc. (pg/mL)	25 400	35 200	105 000	15 140	107 400	21 600	101 800	212 000	210 000	202 000

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		24 Jul 2015