



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07890

Product Description: Cytokine Panel 1 (human) Calibrator Blend
Catalog Number: C0050-2
Lot Number: A0080108
Expiration Date: 30 NOV 2021
Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay									
	hu GM-CSF	hu IL-1 α	hu IL-5	hu IL-7	hu IL-12/23p40	hu IL-15	hu IL-16	hu IL-17A	hu TNF- β	hu VEGF
Reconstituted Calibrator (pg/mL)	1150	468	850	1010	3730	886	3120	6620	693	1340

Reconstitution Volume: 1.0 mL

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. This product 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>	13 FEB 2017