



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07040

Product Description: Cytokine Panel 1 (human) Control Pack

Catalog Number: C4050-1

Lot Number: CP000040

Expiration Date: 30 APR 2020

Concentration of Materials Tested:

Units = pg/mL	Assay									
	hu GM-CSF	hu IL-1 α	hu IL-5	hu IL-7	hu IL-12/23 p40	hu IL-15	hu IL-16	hu IL-17A	hu TNF- β	hu VEGF
Reconstituted Control 1 (A00C0325)	362	167	371	241	1290	241	1160	2590	196	373
Reconstituted Control 2 (A00C0326)	55.8	57.8	123	42.0	208	36.5	328	327	31.7	127
Reconstituted Control 3 (A00C0395)	9.32	19.2	36.0	7.52	37.6	7.94	110	56.9	5.95	46.1

Reconstitution Volume: 250 μ L

Control concentrations were assigned across three kit lots.

Cytokine Panel 1 (human) Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Control 1	A00C0325	2-8°C	30 APR 2020
Cytokine Panel 1 (human) Control 2	A00C0326	2-8°C	31 MAY 2020
Cytokine Panel 1 (human) Control 3	A00C0395	2-8°C	30 APR 2021

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	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	21 JUN 2016