



January 2022

Re: Cytokine Panel 1 (human) Control Pack (Catalog No. C4050-1) Lot No. CP000119, Revised Concentrations

Dear Valued Customer,

As part of our process for continuous product improvement, we periodically review the performance specifications of V-PLEX® products. As we accumulate data from additional lots, if the larger data set indicates a change to a more accurate value, we update the product accordingly.

We recently reviewed the data for the controls in our Cytokine Panel 1 (human) Control Pack. Based on the larger data set, we have decided to reassign the values of the controls for all of the analytes in this panel. We do not believe that the controls have changed. Rather, we believe that the larger data set provides more accurate values for the controls. The revised (Table 1) and original (Table 2) values for the controls are listed below.

Units = pg/mL	Assay									
	GM-CSF	IL-1a	IL-5	IL-7	IL-12/IL-23p40	IL-15	IL-16	IL-17A	TNF-b	VEGF
Control 1 (A00C0774)	425	175	334	323	1200	300	1220	2670	249	354
Control 2 (A00C0775)	66.3	55.6	107	51.7	179	43.1	329	384	40.0	117
Control 3 (A00C0776)	9.52	15.5	32.9	7.03	26.8	7.17	100	52.4	5.35	37.5

Table 1. Revised concentrations

Units = pg/mL	Assay									
	GM-CSF	IL-1a	IL-5	IL-7	IL-12/IL-23p40	IL-15	IL-16	IL-17A	TNF-b	VEGF
Control 1 (A00C0774)	447	173	334	348	1240	314	1160	2650	250	373
Control 2 (A00C0775)	70.5	56.1	108	55.8	189	45.6	327	371	41.6	126
Control 3 (A00C0776)	10.7	16	34.5	7.75	29.9	7.69	100	51.2	5.63	40.1

Table 2. Original concentrations

If you have any questions or concerns, please contact MSD Scientific Support at +1 240-314-2798 or scientificsupport@mesoscale.com.

Sincerely,

MSD Quality Assurance



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Cytokine Panel 1 (human) Control Pack

Catalog Number: C4050-1

Lot Number: CP000119

Expiration Date: 30 APR 2026

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
GM-CSF	425	66.3	9.52	pg/mL
IL-1 α	175	55.6	15.5	pg/mL
IL-5	334	107	32.9	pg/mL
IL-7	323	51.7	7.03	pg/mL
IL-12/23 p40	1200	179	26.8	pg/mL
IL-15	300	43.1	7.17	pg/mL
IL-16	1220	329	100	pg/mL
IL-17A	2670	384	52.4	pg/mL
TNF- β	249	40.0	5.35	pg/mL
VEGF	354	117	37.5	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Control 1	A00C0774	2-8°C	30 APR 2026
Cytokine Panel 1 (human) Control 2	A00C0775	2-8°C	30 APR 2026
Cytokine Panel 1 (human) Control 3	A00C0776	2-8°C	30 APR 2026

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 JAN 2022