



May 2021

Re: Proinflammatory Panel 1 (human) Control Pack (Catalog No. C4049-1) Lot No. CP000114, 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 Proinflammatory 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									
	hu IFN- γ	hu IL-1 β	hu IL-2	hu IL-4	hu IL-6	hu IL-8	hu IL-10	hu IL-12p70	hu IL-13	hu TNF- α
Control 1 (A00C0743)	580	220	382	93.5	259	240	133	171	165	77.8
Control 2 (A00C0744)	161	52.2	39.8	19.1	47.5	37.9	25.0	32.0	45.1	13.7
Control 3 (A00C0745)	42.5	11.0	3.22	2.53	7.38	6.46	3.66	5.72	12.0	2.15

Table 1. Revised concentrations

Units = pg/mL	Assay									
	hu IFN- γ	hu IL-1 β	hu IL-2	hu IL-4	hu IL-6	hu IL-8	hu IL-10	hu IL-12p70	hu IL-13	hu TNF- α
Control 1 (A00C0743)	606	224	385	96.9	263	245	132	176	170	77.9
Control 2 (A00C0744)	167	53.4	40.6	19.6	47.9	38.9	24.9	32.6	45.7	14.2
Control 3 (A00C0745)	46.7	11.7	3.42	2.69	7.72	6.76	3.72	5.99	12.3	2.23

Table 2. Original concentrations

If you have any other 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: Proinflammatory Panel 1 (human) Control Pack

Catalog Number: C4049-1

Lot Number: CP000114

Expiration Date: 31 DEC 2025

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	580	161	42.5	pg/mL
IL-1 β	220	52.2	11.0	pg/mL
IL-2	382	39.8	3.22	pg/mL
IL-4	93.5	19.1	2.53	pg/mL
IL-6	259	47.5	7.38	pg/mL
IL-8	240	37.9	6.46	pg/mL
IL-10	133	25.0	3.66	pg/mL
IL-12p70	171	32.0	5.72	pg/mL
IL-13	165	45.1	12.0	pg/mL
TNF- α	77.8	13.7	2.15	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0743	2-8°C	31 DEC 2025
Proinflammatory Panel 1 (human) Control 2	A00C0744	2-8°C	31 DEC 2025
Proinflammatory Panel 1 (human) Control 3	A00C0745	2-8°C	31 DEC 2025

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		14 MAY 2021