



December 2022

Re: Proinflammatory Panel 1 (human) Control Pack (Catalog No. C4049-1), Lot No. CP000136, 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 Lot No. CP000136. Based on the larger data set, we 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 (A00C0840)	510	223	403	105	265	251	139	183	163	84.1
Control 2 (A00C0841)	136	53.8	39.5	20.9	48.3	39.2	26.4	32.8	43.3	16.8
Control 3 (A00C0842)	40.3	10.6	3.33	3.02	7.51	6.62	3.90	6.42	11.8	2.03

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 (A00C0840)	526	230	419	109	270	261	147	192	169	89.0
Control 2 (A00C0841)	139	54.9	41.0	21.5	48.8	40.6	27.3	34.0	44.3	17.4
Control 3 (A00C0842)	42.0	11.1	3.52	3.18	7.77	6.90	4.02	6.85	12.0	2.06

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: Proinflammatory Panel 1 (human) Control Pack

Catalog Number: C4049-1

Lot Number: CP000136

Expiration Date: 31 JUL 2027

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	510	136	40.3	pg/mL
IL-1 β	223	53.8	10.6	pg/mL
IL-2	403	39.5	3.33	pg/mL
IL-4	105	20.9	3.02	pg/mL
IL-6	265	48.3	7.51	pg/mL
IL-8	251	39.2	6.62	pg/mL
IL-10	139	26.4	3.90	pg/mL
IL-12p70	183	32.8	6.42	pg/mL
IL-13	163	43.3	11.8	pg/mL
TNF- α	84.1	16.8	2.03	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0840	2-8°C	31 JUL 2027
Proinflammatory Panel 1 (human) Control 2	A00C0841	2-8°C	31 JUL 2027
Proinflammatory Panel 1 (human) Control 3	A00C0842	2-8°C	31 JUL 2027

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		12 JAN 2023