



August 2022

Re: Chemokine Panel 1 (human) Control Pack (Catalog No. C4047-1) Lot No. CP000132, 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 Chemokine Panel 1 (human) Control Pack Lot No. CP000132. 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 Eotaxin	hu MIP-1 $\beta$	hu Eotaxin-3	hu TARC	hu IP-10	hu MIP-1 $\alpha$	hu IL-8	hu MCP-1	hu MDC	hu MCP-4
Control 1 (A00C0811)	1770	968	5420	1640	771	1030	61000	470	4660	568
Control 2 (A00C0812)	632	164	647	274	116	397	21400	75.2	2000	226
Control 3 (A00C0813)	195	28.4	93.4	43.3	17.3	165	8060	11.8	937	75.3

Table 1. Revised concentrations

Units = pg/mL	Assay									
	hu Eotaxin	hu MIP-1 $\beta$	hu Eotaxin-3	hu TARC	hu IP-10	hu MIP-1 $\alpha$	hu IL-8	hu MCP-1	hu MDC	hu MCP-4
Control 1 (A00C0811)	1840	941	4700	1780	680	992	55500	500	4940	610
Control 2 (A00C0812)	640	160	615	265	94.9	395	20800	74.9	2010	234
Control 3 (A00C0813)	196	27.6	88.5	43.8	14.3	167	7260	12.4	927	76

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](mailto:scientificsupport@mesoscale.com).

Sincerely,

MSD Quality Assurance



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Control Pack Description:** Chemokine Panel 1 (human) Control Pack

**Catalog Number:** C4047-1

**Lot Number:** CP000132

**Expiration Date:** 30 NOV 2026

**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1770	632	195	pg/mL
MIP-1 $\beta$	968	164	28.4	pg/mL
Eotaxin-3	5420	647	93.4	pg/mL
TARC	1640	274	43.3	pg/mL
IP-10	771	116	17.3	pg/mL
MIP-1 $\alpha$	1030	397	165	pg/mL
IL-8	61000	21400	8060	pg/mL
MCP-1	470	75.2	11.8	pg/mL
MDC	4660	2000	937	pg/mL
MCP-4	568	226	75.3	pg/mL

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0811	2-8°C	30 NOV 2026
Chemokine Panel 1 (human) Control 2	A00C0812	2-8°C	30 NOV 2026
Chemokine Panel 1 (human) Control 3	A00C0813	2-8°C	30 NOV 2026

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		02 SEP 2022