



September 2021

Re: Chemokine Panel 1 (human) Control Pack (Catalog No. C4047-1) Lot No. CP000118, 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. 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									
	Eotaxin	MIP-1β	Eotaxin-3	TARC	IP-10	MIP-1α	IL-8	MCP-1	MDC	MCP-4
Control 1 (A00C0777)	1760	690	3350	2240	599	925	65000	464	3870	670
Control 2 (A00C0766)	634	122	511	278	104	412	22600	72.7	2060	234
Control 3 (A00C0778)	196	20.6	74.3	42.6	15.4	166	8620	12.1	869	74.6

Table 1. Revised concentrations

Units = pg/mL	Assay									
	Eotaxin	MIP-1β	Eotaxin-3	TARC	IP-10	MIP-1α	IL-8	MCP-1	MDC	MCP-4
Control 1 (A00C0777)	1780	710	3200	2290	605	934	65200	478	3890	682
Control 2 (A00C0766)	626	127	489	280	103	425	22800	73.6	2050	237
Control 3 (A00C0778)	194	21.1	74.7	44.5	15.3	172	8020	12.7	867	73.3

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: Chemokine Panel 1 (human) Control Pack 1

Catalog Number: C4047-1

Lot Number: CP000118

Expiration Date: 28 FEB 2026

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1760	634	196	pg/mL
MIP-1 β	690	122	20.6	pg/mL
Exotaxin-3	3350	511	74.3	pg/mL
TARC	2240	278	42.6	pg/mL
IP-10	599	104	15.4	pg/mL
MIP-1 α	925	412	166	pg/mL
hIL-8	65000	22600	8620	pg/mL
MCP-1	464	72.7	12.1	pg/mL
MDC	3870	2060	869	pg/mL
MCP-4	670	234	74.6	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0777	2-8°C	31 MAR 2026
Chemokine Panel 1 (human) Control 2	A00C0766	2-8°C	28 FEB 2026
Chemokine Panel 1 (human) Control 3	A00C0778	2-8°C	31 MAR 2026

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		17 SEP 2021