



December 2021

Re: TH17 Panel 1 (human) Control Pack (Catalog No. C4085-1) Lot No. CP000103, 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 TH17 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 IL-17A Gen. B	Hu IL-21	hu IL-31	hu IL-27	hu IL-23	hu IL-22	hu MIP-3α
Control 1 (A00C0711)	2710	927	985	17200	5440	538	509
Control 2 (A00C0712)	372	203	145	4120	605	88.4	189
Control 3 (A00C0713)	46.5	47.8	21.3	818	65.9	15.4	66.9

Table 1. Revised concentrations

Units = pg/mL	Assay						
	hu IL-17A Gen. B	Hu IL-21	hu IL-31	hu IL-27	hu IL-23	hu IL-22	hu MIP-3α
Control 1 (A00C0711)	2860	1300	971	17500	5210	529	564
Control 2 (A00C0712)	403	297	144	4200	578	88.1	202
Control 3 (A00C0713)	57.1	84.2	22.3	826	63.6	16.6	72.3

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: TH17 Panel 1 (human) Control Pack

Catalog Number: C4085-1

Lot Number: CP000103

Expiration Date: 31 MAR 2025

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IL-17A Gen. B	2710	372	46.5	pg/mL
IL-21	927	203	47.8	pg/mL
IL-31	985	145	21.3	pg/mL
IL-27	17200	4120	818	pg/mL
IL-23	5440	605	65.9	pg/mL
IL-22	538	88.4	15.4	pg/mL
MIP-3 α	509	189	66.9	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
TH17 Panel 1 (human) Control 1	A00C0711	2-8°C	31 MAR 2025
TH17 Panel 1 (human) Control 2	A00C0712	2-8°C	31 MAR 2025
TH17 Panel 1 (human) Control 3	A00C0713	2-8°C	31 MAR 2025

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		10 JAN 2022