

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Chemokine Panel 1 (human) Gen. B Control Pack
Catalog Number: C4705-1
Lot Number: CP000162
Expiration Date: 31 JUL 2029

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Control 1	A00C0965	2-8°C	31 JUL 2029
Chemokine Panel 1 (human) Gen. B Control 2	A00C0958	2-8°C	31 JUL 2029
Chemokine Panel 1 (human) Gen. B Control 3	A00C0956	2-8°C	30 JUN 2029

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1270	198	28.7	pg/mL
MIP-1 β	494	78.4	14.6	pg/mL
Eotaxin-3	13800	1440	125	pg/mL
TARC/CCL17	607	125	29.1	pg/mL
IP-10	1720	193	18.0	pg/mL
MIP-1 α	518	50.5	4.32	pg/mL
MCP-1	338	74.0	18.9	pg/mL
MDC	3700	570	87.7	pg/mL
MCP-4	333	48.8	7.32	pg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		22 AUG 2024