



CERTIFICATE OF ANALYSIS

**Control Pack Description:** Chemokine Panel 1 (human) Gen. B Control Pack  
**Catalog Number:** C4705-1  
**Lot Number:** CP000168  
**Expiration Date:** 30 SEP 2029

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Control 1	A00C0966	2-8°C	30 SEP 2029
Chemokine Panel 1 (human) Gen. B Control 2	A00C0967	2-8°C	30 SEP 2029
Chemokine Panel 1 (human) Gen. B Control 3	A00C0968	2-8°C	30 SEP 2029

**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1340	201	33.7	pg/mL
MIP-1 $\beta$	503	78.8	14.5	pg/mL
Eotaxin-3	14000	1400	124	pg/mL
TARC/CCL17	641	136	24.5	pg/mL
IP-10	1890	209	19.2	pg/mL
MIP-1 $\alpha$	535	48.8	4.29	pg/mL
MCP-1	345	74.9	18.3	pg/mL
MDC	3750	569	170	pg/mL
MCP-4	326	47.5	7.20	pg/mL

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		07 NOV 2024