



CERTIFICATE OF ANALYSIS

Control Pack Description: Chemokine Panel 1 (human) Gen. B Control Pack
Catalog Number: C4705-1
Lot Number: CP000183
Expiration Date: 30 JUN 2029

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Control 1	A00C1023	2-8°C	31 JUL 2030
Chemokine Panel 1 (human) Gen. B Control 2	A00C1024	2-8°C	31 JUL 2030
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	189	28.7	pg/mL
MIP-1 β	556	80.8	14.6	pg/mL
Eotaxin-3	16700	1550	125	pg/mL
TARC/CCL17	593	117	29.1	pg/mL
IP-10	2050	214	18.0	pg/mL
MIP-1 α	521	47.9	4.32	pg/mL
MCP-1	342	74.6	18.9	pg/mL
MDC	3780	554	87.7	pg/mL
MCP-4	331	44.3	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		17 NOV 2025