



CERTIFICATE OF ANALYSIS

Control Pack Description: Chemokine Panel 1 (human) Gen. B Control Pack
Catalog Number: C4705-1
Lot Number: CP000172
Expiration Date: 31 DEC 2029

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Control 1	A00C0980	2-8°C	31 DEC 2029
Chemokine Panel 1 (human) Gen. B Control 2	A00C0993	2-8°C	31 DEC 2029
Chemokine Panel 1 (human) Gen. B Control 3	A00C0994	2-8°C	31 DEC 2029

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1260	180	27.1	pg/mL
MIP-1β	503	76.7	15.2	pg/mL
Eotaxin-3	15800	1490	145	pg/mL
TARC/CCL17	598	120	24.7	pg/mL
IP-10	1740	177	16.4	pg/mL
MIP-1α	567	50.1	4.63	pg/mL
MCP-1	330	76.1	17.6	pg/mL
MDC	3770	552	84.6	pg/mL
MCP-4	321	44.8	8.09	pg/mL

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

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		24 FEB 2025