



CERTIFICATE OF ANALYSIS

Control Pack Description: Chemokine Panel 1 (human) Gen. B Control Pack
Catalog Number: C4705-1
Lot Number: CP000149
Expiration Date: 31 JUL 2028

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Control 1	A00C0912	2-8°C	31 JUL 2028
Chemokine Panel 1 (human) Gen. B Control 2	A00C0913	2-8°C	31 JUL 2028
Chemokine Panel 1 (human) Gen. B Control 3	A00C0914	2-8°C	31 JUL 2028

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1290	180	29.5	pg/mL
MIP-1 β	505	109	27.6	pg/mL
Eotaxin-3	12500	1160	118	pg/mL
TARC/CCL17	579	89.0	16.4	pg/mL
IP-10	1840	183	20.1	pg/mL
MIP-1 α	542	46.7	4.78	pg/mL
MCP-1	347	49.1	7.86	pg/mL
MDC	3920	527	85.3	pg/mL
MCP-4	339	46.9	7.62	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		27 SEP 2023