

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Chemokine Panel 1 (human) Gen. B Control Pack
Catalog Number: C4705-1
Lot Number: CP000192
Expiration Date: 31 MAR 2031

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Control 1	A00C1056	2-8°C	31 MAR 2031
Chemokine Panel 1 (human) Gen. B Control 2	A00C1057	2-8°C	30 APR 2031
Chemokine Panel 1 (human) Gen. B Control 3	A00C1058	2-8°C	30 APR 2031

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1260	188	32.7	pg/mL
MIP-1 β	540	89.2	19.2	pg/mL
Eotaxin-3	12800	1310	139	pg/mL
TARC/CCL17	624	123	26.6	pg/mL
IP-10	1860	198	23.8	pg/mL
MIP-1 α	531	48.8	5.31	pg/mL
MCP-1	331	77.4	20.7	pg/mL
MDC	3680	563	102	pg/mL
MCP-4	317	44.5	8.19	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		26 JUN 2026