



CERTIFICATE OF ANALYSIS

Control Pack Description: Chemokine Panel 1 (human) Control Pack
Catalog Number: C4047-1
Lot Number: CP000178
Expiration Date: 30 SEP 2029

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0990	2-8°C	31 DEC 2029
Chemokine Panel 1 (human) Control 2	A00C1004	2-8°C	31 MAR 2030
Chemokine Panel 1 (human) Control 3	A00C0964	2-8°C	30 SEP 2029

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1740	622	167	pg/mL
MIP-1 β	761	128	28.4	pg/mL
Eotaxin-3	4290	612	90.1	pg/mL
TARC/CCL17	1320	219	39.3	pg/mL
IP-10	637	100	20.2	pg/mL
MIP-1 α	966	416	177	pg/mL
IL-8 (HA)	60200	21500	7770	pg/mL
MCP-1	467	68.1	12.5	pg/mL
MDC	9810	1750	1880	pg/mL
MCP-4	769	207	62.9	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		03 JUN 2025