



CERTIFICATE OF ANALYSIS

Control Pack Description: Chemokine Panel 1 (human) Control Pack
Catalog Number: C4047-1
Lot Number: CP000171
Expiration Date: 30 SEP 2029

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0992	2-8°C	31 DEC 2029
Chemokine Panel 1 (human) Control 2	A00C0991	2-8°C	31 DEC 2029
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	1670	583	167	pg/mL
MIP-1 β	741	125	28.4	pg/mL
Eotaxin-3	3780	620	90.1	pg/mL
TARC/CCL17	1070	199	39.3	pg/mL
IP-10	565	94.0	20.2	pg/mL
MIP-1 α	925	407	177	pg/mL
IL-8 (HA)	57400	19500	7770	pg/mL
MCP-1	428	67.6	12.5	pg/mL
MDC	8770	4070	1880	pg/mL
MCP-4	737	230	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		07 FEB 2025