



CERTIFICATE OF ANALYSIS

Control Pack Description: Chemokine Panel 1 (human) Control Pack
Catalog Number: C4047-1
Lot Number: CP000152
Expiration Date: 30 SEP 2028

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0921	2-8°C	30 SEP 2028
Chemokine Panel 1 (human) Control 2	A00C0922	2-8°C	30 SEP 2028
Chemokine Panel 1 (human) Control 3	A00C0923	2-8°C	30 SEP 2028

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1720	577	176	pg/mL
MIP-1 β	773	136	24.1	pg/mL
Eotaxin-3	4400	648	110	pg/mL
TARC/CCL17	1450	230	37.7	pg/mL
IP-10	737	116	19.9	pg/mL
MIP-1 α	995	441	172	pg/mL
IL-8 (HA)	64700	21700	8320	pg/mL
MCP-1	475	72.2	12.1	pg/mL
MDC	4700	2060	949	pg/mL
MCP-4	655	207	63.7	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		30 JUL 2024