



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Chemokine Panel 1 (human) Control Pack

Catalog Number: C4047-1

Lot Number: CP000107

Expiration Date: 31 JUL 2025

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1770	650	204	pg/mL
MIP-1 β	1100	164	26.5	pg/mL
Eotaxin-3	4000	577	89.0	pg/mL
TARC	1440	235	39.8	pg/mL
IP-10	580	85.6	14.5	pg/mL
MIP-1 α	931	402	168	pg/mL
IL-8	63800	22600	8010	pg/mL
MCP-1	476	70.9	13.6	pg/mL
MDC	8920	2750	892	pg/mL
MCP-4	671	248	76.2	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0726	2-8°C	31 JUL 2025
Chemokine Panel 1 (human) Control 2	A00C0727	2-8°C	31 JUL 2025
Chemokine Panel 1 (human) Control 3	A00C0728	2-8°C	31 JUL 2025

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 OCT 2020