



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4047-1

Lot Number: CP000082

Expiration Date: 30 NOV 2023

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1610	590	182	pg/mL
MIP-1 β	1040	178	27.1	pg/mL
Eotaxin-3	4300	615	95.7	pg/mL
TARC	1340	223	36.9	pg/mL
IP-10	561	85.6	14	pg/mL
MIP-1 α	911	401	166	pg/mL
IL-8	61100	21300	8370	pg/mL
MCP-1	465	71.3	11.5	pg/mL
MDC	8960	2630	645	pg/mL
MCP-4	660	239	71.2	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0618	2-8°C	30 NOV 2023
Chemokine Panel 1 (human) Control 2	A00C0619	2-8°C	30 NOV 2023
Chemokine Panel 1 (human) Control 3	A00C0620	2-8°C	30 NOV 2023

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		10 SEP 2020