



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C4047-1

**Lot Number:** CP000076

**Expiration Date:** 30 JUN 2023


### Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1550	594	182	pg/mL
MIP-1 $\beta$	991	172	25.5	pg/mL
Eotaxin-3	4400	654	99.6	pg/mL
TARC	1300	216	35.4	pg/mL
IP-10	544	86.4	14.2	pg/mL
MIP-1 $\alpha$	926	404	162	pg/mL
IL-8(HA)	63 500	21 900	8970	pg/mL
MCP-1	465	71.7	12.1	pg/mL
MDC	8970	2710	907	pg/mL
MCP-4	650	246	79.2	pg/mL

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0588	2-8°C	30 JUN 2023
Chemokine Panel 1 (human) Control 2	A00C0589	2-8°C	30 JUN 2023
Chemokine Panel 1 (human) Control 3	A00C0590	2-8°C	30 JUN 2023

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality		30 APR 2019