



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Chemokine Panel 1 (human) Control Pack

Catalog Number: C4047-1

Lot Number: CP000064

Expiration Date: 31 OCT 2022

Concentration of Materials Tested:

Units = pg/mL	Assay									
	hu Eotaxin	hu MIP-1 β	hu Eotaxin-3	hu TARC	hu IP-10	hu MIP-1 α	hu IL-8	hu MCP-1	hu MDC	hu MCP-4
Control 1 (A00C0525)	1640	734	3705	1280	364	824	139500	568	11000	762
Control 2 (A00C0526)	558	116	526	184	55	330	60750	83	2965	280
Control 3 (A00C0527)	170.5	17.1	74.0	30.8	9.4	125.0	24500	13.8	869.0	87.3

Reconstitution Volume: 250 μ L

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0525	2-8°C	31 OCT 2022
Chemokine Panel 1 (human) Control 2	A00C0526	2-8°C	31 OCT 2022
Chemokine Panel 1 (human) Control 3	A00C0527	2-8°C	31 OCT 2022

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	19 DEC 2017