



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Chemokine Panel 1 (human) Control Pack

Catalog Number: C4047-1

Lot Number: CP000053

Expiration Date: 30 NOV 2021

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 (A00C0442)	1410	904	4870	1360	473	795	128000	407	8530	614
Control 2 (A00C0443)	536	156	816	232	74.9	362	59400	65.9	2680	234
Control 3 (A00C0444)	177	24.5	137	38.3	11.7	140	23700	11.6	786	71.2

Reconstitution Volume: 250 μ L

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0442	2-8°C	30 NOV 2021
Chemokine Panel 1 (human) Control 2	A00C0443	2-8°C	30 NOV 2021
Chemokine Panel 1 (human) Control 3	A00C0444	2-8°C	30 NOV 2021

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>	04 MAY 2017