



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Chemokine Panel 1 (human) Control Pack

Catalog Number: C4047-1

Lot Number: CP000066

Expiration Date: 31 JAN 2023

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 (A00C0547)	1580	940	3920	1390	574	887	70 100	595	10 800	688
Control 2 (A00C0548)	311	158	553	213	87.5	393	23 500	155	3120	258
Control 3 (A00C0566)	171	26.9	86.3	37.8	15.4	152	9330	15.2	982	78.2

Reconstitution Volume: 250 μ L

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0547	2-8°C	31 JAN 2023
Chemokine Panel 1 (human) Control 2	A00C0548	2-8°C	31 JAN 2023
Chemokine Panel 1 (human) Control 3	A00C0566	2-8°C	31 JAN 2023

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>	23 FEB 2018