



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Chemokine Panel 1 (human) Control Pack

Catalog Number: C4047-1

Lot Number: CP000058

Expiration Date: 29 FEB 2020

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 (A00C0498)	1370	797	4760	1130	348	824	171 000	420	11 100	635
Control 2 (A00C0499)	518	130	755	197	57.3	371	58 100	66.6	3460	232
Control 3 (A00C0500)	160	22.4	118	34.2	10.3	152	22 500	11.1	907	82.2

Reconstitution Volume: 250 μ L

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0498	2-8°C	30 APR 2022
Chemokine Panel 1 (human) Control 2	A00C0499	2-8°C	29 FEB 2020
Chemokine Panel 1 (human) Control 3	A00C0500	2-8°C	30 APR 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>	08 AUG 2017