



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05656

Product Description: Chemokine Panel 1 (human) Control Pack

Catalog Number: C4047-1

Lot Number: CP000028

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
Reconstituted Control 1 (A00C0304)	1230	768	4670	1160	381	826	150000	437	8670	601
Reconstituted Control 2 (A00C0305)	443	124	794	188	61.8	369	53200	71.5	2660	216
Reconstituted Control 3 (A00C0306)	140	22.4	127	34.3	10.8	149	20400	11.6	703	75.0

Reconstitution Volume: 250 μ L

Control concentrations were assigned across three kit lots.

Chemokine Panel 1 (human) Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C0304	2-8°C	29 FEB 2020
Chemokine Panel 1 (human) Control 2	A00C0305	2-8°C	29 FEB 2020
Chemokine Panel 1 (human) Control 3	A00C0306	2-8°C	31 MAR 2020

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Michael Grossman	Quality Assurance		25 JUN 2015