



# CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-04712

**Product Description:** Chemokine Panel 1 (human) Control Pack

**Catalog Number:** C4047-1

**Lot Number:** CP000021

**Expiration Date:** 31 JUL 2019

## Concentration of Materials Tested:

Units = pg/mL	Assay									
	hu Eotaxin	hu MIP-1 $\beta$	hu Eotaxin-3	hu TARC	hu IP-10	hu MIP-1 $\alpha$	hu IL-8	hu MCP-1	hu MDC	hu MCP-4
Reconstituted Control 1 (A00C0267)	1040	603	3720	971	1470	709	161000	382	5790	475
Reconstituted Control 2 (A00C0268)	283	156	897	235	378	188	52300	88.9	1330	131
Reconstituted Control 3 (A00C0269)	65.6	36.7	226	63.4	93.5	42.8	16100	22.7	333	32.2

**Reconstitution Volume:** 250  $\mu$ 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	A00C0267	2-8 $^{\circ}$ C	31 JUL 2019
Chemokine Panel 1 (human) Control 2	A00C0268	2-8 $^{\circ}$ C	31 JUL 2019
Chemokine Panel 1 (human) Control 3	A00C0269	2-8 $^{\circ}$ C	31 JUL 2019

**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
<b>Review/Approval</b>	Caitlin Hansen	Quality Control		27 AUG 2014