



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05081

Product Description: Chemokine Panel 1 (human) Calibrator Blend

Catalog Number: C0047-2

Lot Number: A0080046

Expiration Date: 31 MAY 2017

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	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 Calibrator (pg/mL)	1440	982	4760	1550	2470	1010	68 100	445	10 100	623

Reconstitution Volume: 1.0 mL

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	Caitlin Hansen	Quality Control		