



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Chemokine Panel 1 (human) Calibrator Blend

Catalog Number: C0047-2

Lot Number: A0080183

Expiration Date: 31 OCT 2023

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)	1960	1240	5960	1800	2480	1090	72 100	537	10 200	751

Reconstitution Volume: 1000 μ L

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>	16 NOV 2018