



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Chemokine Panel 1 (human) Calibrator Blend

**Catalog Number:** C0047-2

**Lot Number:** A0080131

**Expiration Date:** 31 JUL 2022

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

	Assay									
	hu Eotaxin	hu MIP-1 $\beta$	hu Exotaxin-3	hu TARC	hu IP-10	hu MIP-1 $\alpha$	hu IL-8	hu MCP-1	hu MDC	hu MCP-4
Reconstituted Calibrator (pg/mL)	1860	1050	5410	1630	2420	1030	86 200	686	12 200	787

**Reconstitution Volume:** 1000  $\mu$ L

**Statement:** This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	31 JUL 2017