



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0047-2

**Lot Number:** A0080122

**Expiration Date:** 28 FEB 2022

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

	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 Calibrator (pg/mL)	1840	1110	7120	1890	2660	1080	85 700	517	10 800	728

**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>	12 JAN 2018