



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Cytokine Panel 1 (mouse) Calibrator Blend

**Catalog Number:** C0245-2

**Lot Number:** A0080218

**Expiration Date:** 31 OCT 2024

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-9	3720	pg/mL
MCP-1	447	pg/mL
IL-33	3010	pg/mL
IL-27p28/IL-30	8830	pg/mL
IL-15	39400	pg/mL
IL-17A/F	2380	pg/mL
MIP-1 $\alpha$	739	pg/mL
IP-10	1100	pg/mL
MIP-2	559	pg/mL

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

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Carissa Hubbard	Quality		26 OCT 2019