



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0245-2

**Lot Number:** A0080156

**Expiration Date:** 31 MAR 2023

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

	Assay								
	ms IL-9	ms MCP-1	ms IL-33	ms IL-27p28/ IL-30	ms IL-15	ms IL17A/F	ms MIP-1 $\alpha$	ms IP-10	ms MIP-2
Reconstituted Calibrator (pg/mL)	3730	426	3050	7940	36 000	2370	601	896	596

**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>	01 JUN 2018