



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0245-2

**Lot Number:** A0080135

**Expiration Date:** 30 JUN 2022

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

	Assay								
	ms IL-9	ms MCP-1	ms IL-33	ms IL-27p78/ IL-30	ms IL-15	ms IL-17A/F	ms MIP-1 $\alpha$	ms IP-10	ms MIP-2
Reconstituted Calibrator (pg/mL)	3970	482	3320	10400	38100	2530	619	910	635

**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>	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	26JUN2017