



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 1 (mouse) Calibrator Blend

Catalog Number: C0245-2

Lot Number: A0080125

Expiration Date: 31 MAR 2022

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 IL-17A/F	ms MIP-1a	ms IP-10	ms MIP-2
Reconstituted Calibrator (pg/mL)	3860	491	3280	10 500	37 600	2460	601	898	646

Reconstitution Volume: 1000 µL

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	06 JUN 2017