



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0050-2

**Lot Number:** A0080188

**Expiration Date:** 30 NOV 2023

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

	Assay									
	hu GM-CSF	hu IL-1 $\alpha$	hu IL-5	hu IL-7	hu IL-12/23p40	hu IL-15	hu IL-16	hu IL-17A	hu TNF- $\beta$	hu VEGF
Reconstituted Calibrator (pg/mL)	1145	454	868	859	3485	877	3075	5740	646	1080

**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>	20 DEC 2018