



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Cytokine Panel 2 (human) Calibrator Blend

**Catalog Number:** C0084-2

**Lot Number:** A0080116

**Expiration Date:** 31 JAN 2022

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

	Assay							
	hu IL-17A/F	hu IL-17B	hu IL-17C	hu IL-17D	hu IL-1RA	hu IL-3	hu IL-9	hu TSLP
Reconstituted Calibrator (pg/mL)	6140	1400	2500	7240	714	3430	1440	515

**Reconstitution Volume:** 1.0 mL

**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>	21 JUL 2017