



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 2 (human) Calibrator Blend

Catalog Number: C0084-2

Lot Number: A0080126

Expiration Date: 31 MAR 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)	5607	1520	2627	7467	732	3060	1490	563

Reconstitution Volume: 1.00 mL

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

	Name	Function	Signature	Date
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	13JUN2017