



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Proinflammatory Panel 1 (rat) Calibrator Blend

**Catalog Number:** C0044-2

**Lot Number:** A0080124

**Expiration Date:** 31 MAR 2022

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

	Assay								
	rt IFN- $\gamma$	rt IL-1 $\beta$	rt IL-4	rt IL-5	rt IL-6	rt KC/GRO	rt IL-10	rt IL-13	rt TNF- $\alpha$
Reconstituted Calibrator (pg/mL)	6810	12 800	1310	10 500	14 600	3360	25 000	1780	1360

**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>	18 SEP 2017