



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0048-2

**Lot Number:** A0080197

**Expiration Date:** 31 DEC 2023

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

	Assay									
	ms IFN- $\gamma$	ms IL-1 $\beta$	ms IL-2	ms IL-4	ms IL-5	ms IL-6	ms KC/GRO	ms IL-10	ms IL-12p70	ms TNF- $\alpha$
Reconstitution Calibrator (pg/mL)	949	1 480	2 450	1 570	892	4 660	2 160	2 510	31 900	577

**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>	24 JAN 2019