



### Certificate of Analysis

#### VeriKine-HS™ Pig Interferon Alpha ELISA Kit

Catalog No: 47100-1

Lot No: 7177

Expiration: January 31, 2021

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP256	K5894	1
Plate Sealers	N/A	N/A	4
Wash Solution Concentrate	SMP022-60	K5832	50 ml
Pig IFN Alpha Standard 10,000 pg/ml	SMP257-1	K5897	1 vial
Dilution Buffer	SMP021-15	K5511	15 ml
Assay Buffer	SMP258-15	K5898	15 ml
Antibody Concentrate	SMP259-1	K5899	1 vial
Antibody Diluent	SMP260-15	K5901	15 ml
HRP Conjugate Concentrate	SMP056-240	K5900	1 vial
Concentrate Diluent	SMP024-15	K5743	15 ml
TMB Substrate Solution	KET-15	180803D02	15 ml
Stop Solution	SCY-15	51066	15 ml

#### Lot Specific Preparation

(Refer to VeriKine-HS™ Pig IFN α ELISA Kit Protocol for full instructions.)

#### Antibody Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
<b>Antibody Concentrate (µl)</b>	20	40	60	80	100	120
<b>Antibody Diluent (ml)</b>	2	4	6	8	10	12

#### HRP Solution Preparation

Micro-plate Strips Used	2	4	6	8	10	12
<b>HRP Conjugate Concentrate (µl)</b>	20	40	60	80	100	120
<b>Concentrate Diluent (ml)</b>	2	4	6	8	10	12

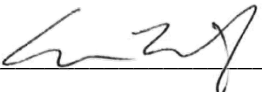
#### Performance Characteristics

Intra-Assay CV: ≤ 8%

Inter-Assay CV: ≤ 10%

Average Spike Recovery: 107%

#### Authorization

Released by: 

Date: February 4, 2020

Sold under license from Pestka Biomedical Laboratory, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to humans. Not for resale in original or any modified form, including inclusion in a kit for any purpose. Not for use in the preparation of any commercial product.