



### Certificate of Analysis

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

Catalog No: 47100-1

Lot No: 7231

Expiration: May 31, 2021

Kit Components	Part No.	Lot No.	Quantity
Plate(s)	SMP256	K6096	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	K6087	1 vial
Dilution Buffer	SMP021-15	K5151	15 ml
Assay Buffer	SMP258-15	K6088	15 ml
Antibody Concentrate	SMP259-1	K6089	1 vial
Antibody Diluent	SMP260-15	K6091	15 ml
HRP Conjugate Concentrate	SMP056-240	K6095	1 vial
Concentrate Diluent	SMP024-15	K5743	15 ml
TMB Substrate Solution	KET-15	180803D02	15 ml
Stop Solution	SCY-15	56334	15 ml

#### Lot Specific Preparation

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

#### Antibody Solution Preparation

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

#### Performance Characteristics

Intra-Assay CV:  $\leq$  8%

Inter-Assay CV:  $\leq$  10%

Average Spike Recovery: 107%

#### Authorization

Released by: \_\_\_\_\_

Date: May 25, 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.