

Research Use Only. Not for use in diagnostic procedures.

PD-1 and PD-L1 (Cynomolgus) Binding AlphaLISA Kit

Product No.: AL583 C/F

Contents

	Page
Product Information.....	2
Quality Control.....	2
Analyte of Interest.....	3
Description of the AlphaLISA Assay	3
Precautions.....	3
Kit content: Reagents and Materials.....	4
Recommendations.....	5
Competition Assay Procedure.....	5
Troubleshooting Guide.....	10

Product Information

Application: This kit is designed to assess inhibitors of cynomolgus PD-1 and cynomolgus PD-L1 binding, using a homogeneous AlphaLISA assay (no wash steps). This assay can facilitate the design and development of antibody therapeutics by using competitive binding to cynomolgus PD-1/PD-L1 to complement PerkinElmer's PD-1/PD-L1 (Human) AlphaLISA Binding kit (Cat# AL356) and PD-1/PD-L1 (Mouse) AlphaLISA Binding Kit (Cat# AL580)

Sensitivity: IC_{50} : 0.062 μ g/mL (average, using anti hPD-L1 antibody, BioLegend Cat # 329926)

Signal to background ratio: 910 using 5 nM cynoPD-1 and 5 nM cynoPD-L1

Kit contents: The kit contains 5 components: anti-6xHis AlphaLISA Acceptor beads, Streptavidin-coated Donor beads, Biotinylated cynomolgus PD-L1, His tagged cynomolgus PD-1 and 10X Binding Assay Buffer.

Storage: The kit components must be stored at 4 °C in the dark. Reconstituted proteins can be stored at –20 °C for 3 months.

Stability: This kit is stable for at least 6 months from the manufacturing date when stored in its original packaging and the recommended storage conditions.

Quality Control

Lot to lot consistency is confirmed in an AlphaLISA assay. Maximum and minimum signals were measured on the EnVision Multilabel Plate Reader with Alpha option using the protocol described in this technical data sheet. We certify that these results meet our quality release criteria. Maximum and minimum counts may vary between bead lots and the instrument used.

Analyte of Interest

Programmed cell death protein 1 (PD-1), also known as cluster of differentiation 279 (CD279), belongs to immunoglobulin superfamily and is a transmembrane receptor protein. Programmed death ligand 1 (PD-L1), also known as cluster of differentiation 274 (CD274) or B7 homolog1 (B7-H1) belongs to the growing B7 family of immune proteins. Cynomolgus PD-1 shares 95% with human PD-1 extracellular domain (ECD) and cynomolgus PD-L1 shares 92% with human PD-L1 ECD. PD-L1 and PD-L2 are two ligands for PD-1. By binding to PD-1 on activated T-cells and B-cells, PD-L1 may inhibit ongoing T-cell responses by inducing apoptosis and arresting cell-cycle progression. Accordingly, it leads to growth of immunogenic tumors by increasing apoptosis of antigen specific T cells and may contribute to immune evasion by cancers. Therefore, blocking PD-1 and PD-L1 binding is considered as a promising therapeutic target for autoimmune disease and malignant cancers.

Description of the AlphaLISA Assay

The AlphaLISA detection of cynomolgus (cyno) PD-1/PD-L1 binding uses anti-6xHis AlphaLISA® acceptor beads to capture the His tagged cynoPD-1 and Streptavidin-coated donor beads to capture the biotinylated cynoPD-L1. Donor beads and acceptor beads come into proximity through cynoPD-L1 binding to cynoPD-1. Excitation of the Donor beads provokes the release of singlet oxygen that triggers a cascade of energy transfer reactions in the Acceptor beads, resulting in a sharp emission at 615 nm (Figure 1).

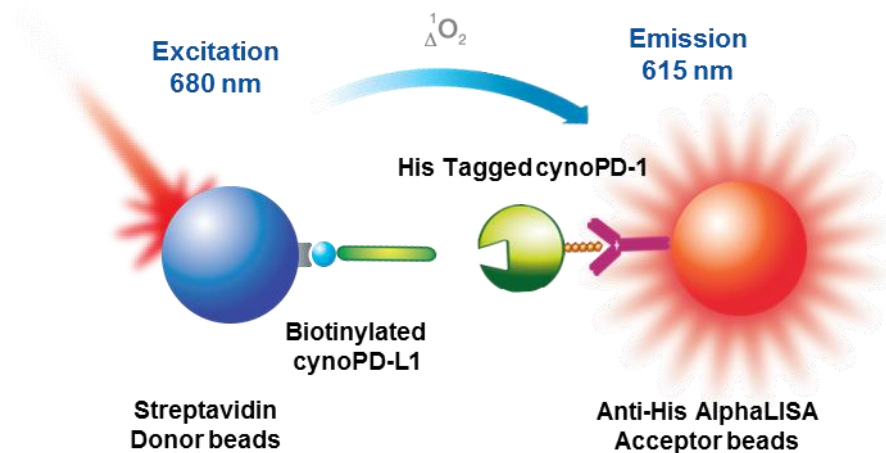


Figure 1. AlphaLISA Assay Principle.

Precautions

- The Alpha Donor beads are light-sensitive. All the other assay reagents can be used under normal light conditions. All Alpha assays using the Donor beads should be performed under subdued laboratory lighting (< 100 lux). Green filters (LEE 090 filters (preferred) or Roscolux filters #389 from Rosco) can be applied to light fixtures.
- All blood components and biological materials should be handled as potentially hazardous. The proteins included in this kit are from a human source.

Kit Content: Reagents and Materials

Kit components	AL583C*** (500 assay points)	AL583F*** (5000 assay points)
Anti-6xHis AlphaLISA Acceptor beads stored in PBS, 0.05% Kathon CG/ICP, pH 7.2	20 µL @ 5 mg/mL (1 brown tube, <u>white</u> cap)	200 µL @ 5 mg/mL (1 brown tube, <u>white</u> cap)
Streptavidin (SA)-coated Donor beads stored in 25 mM HEPES, 100 mM NaCl, 0.05% Kathon CG/ICP, pH 7.4	20 µL @ 5 mg/mL (1 brown tube, <u>black</u> cap)	200 µL @ 5 mg/mL (1 brown tube, <u>black</u> cap)
Lyophilized cynoPD-L1 (Biotinylated)*	2.59 µg, lyophilized (1 tube, <u>clear</u> cap)	2.59 µg, lyophilized (10 tubes, <u>clear</u> caps)
Lyophilized cynoPD-1 (His tagged)*	0.84 µg, lyophilized (1 tube, <u>clear</u> cap)	0.84 µg, lyophilized (10 tubes, <u>clear</u> caps)
AlphaLISA 10X Binding Assay Buffer**	10 mL, 1 small bottle	100 mL, 1 large bottle

* Reconstitute cynoPD-1 and cynoPD-L1 in 100 µL Milli-Q® grade H₂O respectively. The reconstituted proteins should be used within 60 minutes or aliquoted into screw-capped polypropylene vials and stored at -20 °C for further experiments. Avoid multiple freeze-thaw cycles.

** Extra buffer can be ordered separately (cat # AL018C: 10 mL, cat # AL018F: 100 mL).

*** The number of assay points is based on an assay volume of 20 µL in 384 well plates using the kit components at the recommended concentrations (5 nM).

Sodium azide should **not** be added to the stock reagents. High concentrations of sodium azide (> 0.001 % final in the assay) might decrease the AlphaLISA signal.

Specific additional required reagents and materials:

The following materials are recommended:

Item	Suggested source	Catalog #
TopSeal™-A Adhesive Sealing Film	PerkinElmer Inc.	6050185
AlphaPlate-384, Shallow Well (ProxiPlate)	PerkinElmer Inc.	6008350 6008359
EnVision®-Alpha Reader	PerkinElmer Inc.	-

The following reagents might be required for particular applications:

Item	Supplier	Catalog number
Anti-human PD-1 antibody	BioLegend	329926
Mouse IgG1 k, control	BioLegend	400165
Anti-human PD-L1 antibody	R&D Systems	MAB1562
Rabbit IgG, control	Jackson Immuno Research	011-000-003
Cyno PD-1	AcroBiosystems	PD1-C5254
Cyno PD-L1	AcroBiosystems	PD1-C5253

Recommendations

- The volume indicated on each tube is guaranteed for single pipetting. Multiple pipetting of the reagents may reduce the theoretical amount left in the tube. To minimize loss when pipetting beads, it is preferable not to pre-wet the tip.
- Centrifuge all tubes (including lyophilized proteins) before use to improve recovery of content (2000g, 10-15 sec). Re-suspend the beads by vortexing before use. Do not vortex the proteins.
- Use Milli-Q® grade H₂O to reconstitute the lyophilized proteins.
- When reagents are added to the microplate, make sure the liquids are at the bottom of the well.
- Small volume may be prone to evaporation. It is recommended to cover microplates with TopSeal™-A Plus Adhesive Sealing Films to reduce evaporation during incubation. Microplates can be read with the TopSeal™-A Film.
- The AlphaLISA signal is detected with an EnVision Multilabel Reader equipped with the Alpha option using the AlphaScreen standard settings (e.g. Total Measurement Time: 550 ms, Laser 680 nm Excitation Time: 180 ms, Mirror: D640as, Emission Filter: M570w, Center Wavelength 570 nm, Bandwidth 100 nm, Transmittance 75%).
- AlphaLISA signal will vary with temperature and incubation time. For consistent results, identical incubation times and temperature should be used for each plate.

Competition Assay Procedure

IMPORTANT: PLEASE READ THE RECOMMENDATIONS BELOW BEFORE USE

- The protocol described below is an **example** for generating 500 assay points in a 20 µL final assay volume per point. If a different amount of samples are tested, the volumes of all reagents have to be adjusted accordingly. These calculations do not include excess reagent to account for losses during transfer of solutions or dead volumes.
- The dilution protocol is provided for information only. As needed, the number of replicates or the range of concentrations covered can be modified.

One Incubation Step Protocol described as below:

1) Preparation of 1X Binding Assay Buffer:

Add 1 mL of 10X Binding Assay Buffer to 9 mL Milli-Q® grade H₂O.

2) Serial dilutions of 4X anti PD-1 or anti PD-L1 antibody in 1X Binding Assay Buffer as follows:

Tube	Volume of Antibody	Volume of buffer	[Ab] (g/mL) (4X)	[Ab] (g/mL) (1X)
A	4 µL of 1 mg/mL stock	96 µL	4.00E-05	1.00E-05
B	30 µL of tube A	70 µL	1.20E-05	3.00E-06
C	30 µL of tube B	60 µL	4.00E-06	1.00E-06
D	30 µL of tube C	70 µL	1.20E-06	3.00E-07
E	30 µL of tube D	60 µL	4.00E-07	1.00E-07
F	30 µL of tube E	70 µL	1.20E-07	3.00E-08
G	30 µL of tube F	60 µL	4.00E-08	1.00E-08
H	30 µL of tube G	70 µL	1.20E-08	3.00E-09
I	30 µL of tube H	60 µL	4.00E-09	1.00E-09
J	30 µL of tube I	70 µL	1.20E-09	3.00E-10
K	30 µL of tube J	60 µL	4.00E-10	1.00E-10
L	0	70 µL	0	0

3) Preparation of 4X his tagged cynoPD-1 (20 nM):

- Reconstitute lyophilized cynoPD-1 (0.84 µg) in 100 µL H₂O to make 500 nM cynoPD-1.
- Add 100 µL of 500 nM cynoPD-1 to 2400 µL 1X Binding Assay Buffer.
- Prepare just before use and diluted cynoPD-1 should be added to the assay plate in 10 minutes.

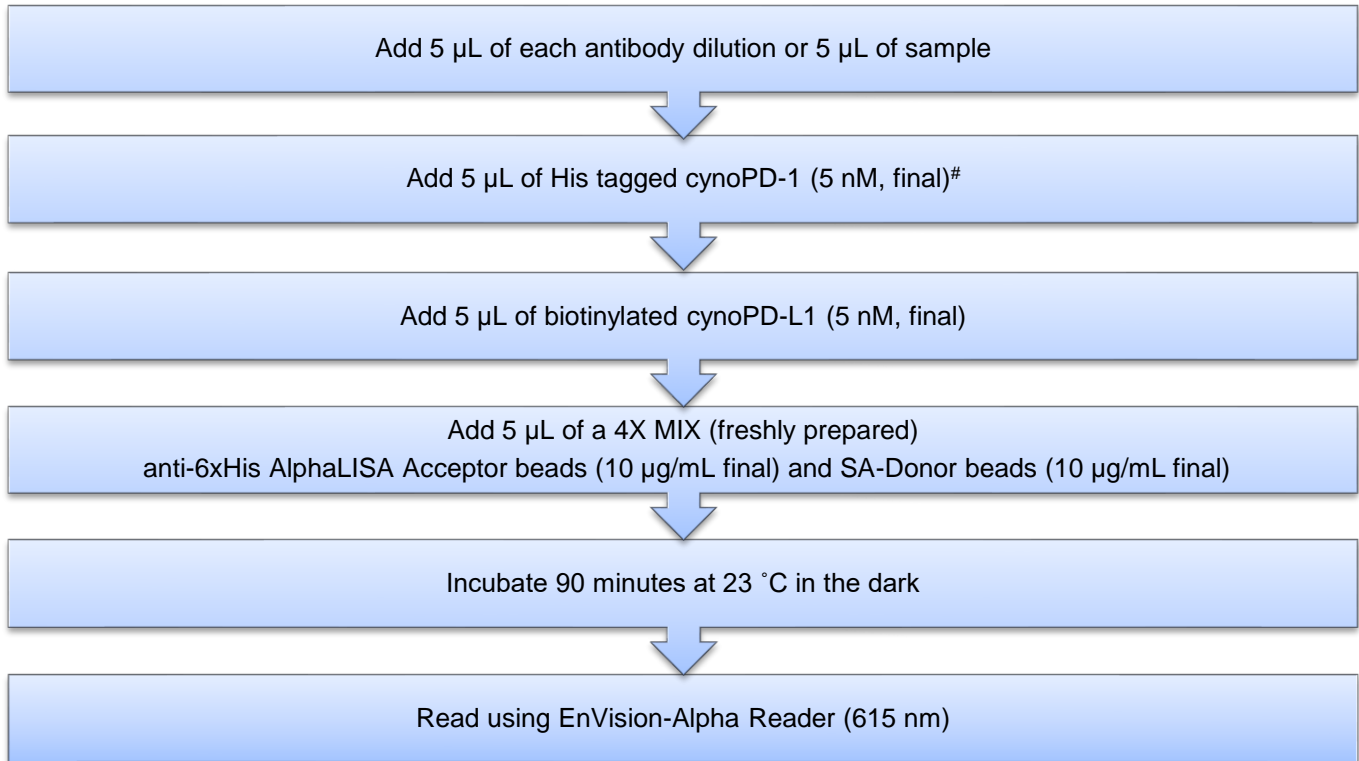
4) Preparation of 4X biotinylated cynoPD-L1 (20 nM):

- Reconstitute lyophilized cynoPD-L1 (2.59 µg) in 100 µL H₂O to make 500 nM cynoPD-L1.
- Add 100 µL of 500 nM cynoPD-L1 to 2400 µL 1X Binding Assay Buffer.
- Prepare just before use and diluted cynoPD-L1 should be added to the assay plate in 10 minutes.

5) Preparation of 4X mix of Anti-6xHis AlphaLISA Acceptor beads (40 µg/mL) and Streptavidin (SA) Donor beads (40 µg/mL):

- Keep the beads under subdued laboratory lighting.
- Add 20 µL of 5 mg/mL Anti-6xHis AlphaLISA Acceptor beads and 20 µL of 5 mg/mL SA-Donor beads to 2460 µL of 1X Binding Assay Buffer.
- Prepare just before use.

6) In a Shallow ProxiPlate (384 wells):



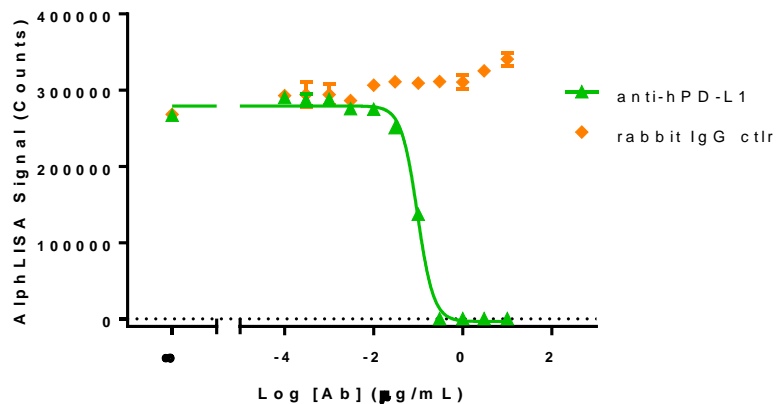
Read Settings: AlphaLISA signal is detected using an EnVision Multilabel Reader equipped with the Alpha option using the following settings: Total Measurement Time: 550 ms, Laser: 680 nm, Excitation Time: 180 ms, Mirror: 640as (Barcode# 444), Emission Filter: Wavelength 570nm, bandwidth: 100nm, Transmittance 75%, (Barcode# 244).

If screening anti-cynoPD-L1 antibodies, add cynoPD-L1 first, then add cynoPD-1.

Typical competitive binding Data:

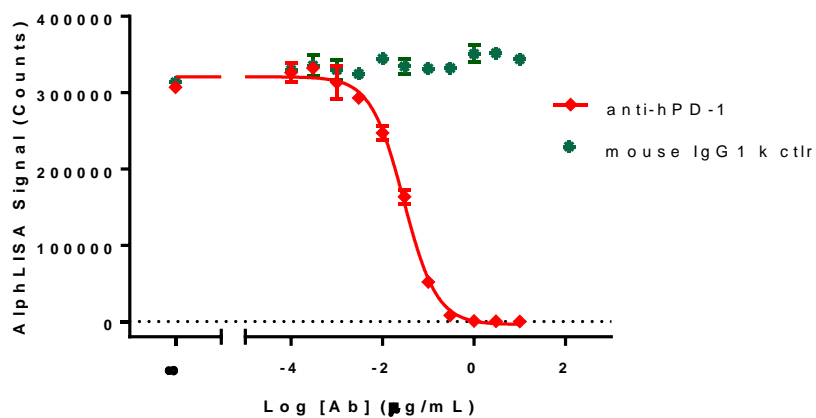
(A)

Anti-hPD-L1 antibody blocking binding to cynoPD-1:PD-L1 (5nM:5nM)



(B)

Anti-hPD-1 antibody blocking binding to cynoPD-1:PD-L1 (5nM:5nM)



(C)

Unlabeled protein competitive binding to cynoPD-1:PD-L1 (5 nM :5 nM)

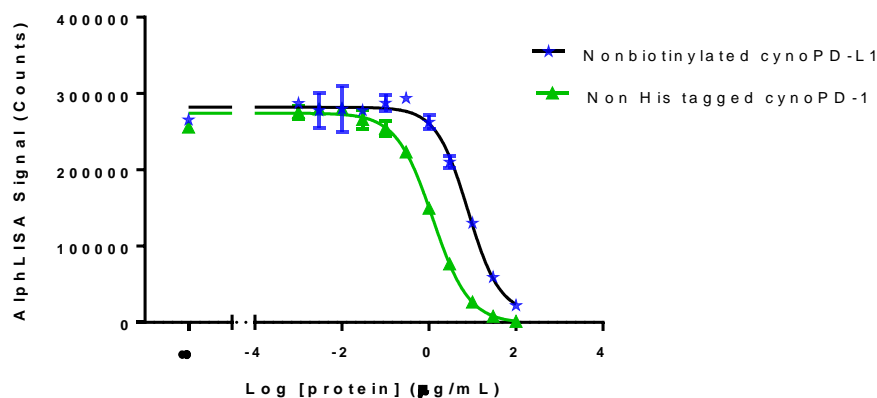


Figure 2. Competitive Binding: (A) Anti-hPD-L1 antibody blocking cynoPD-1/PD-L1 binding with $IC_{50} = 0.096 \mu\text{g/mL}$. Rabbit IgG was measured as a negative control. (B) Anti-hPD-1 antibody blocking cynoPD-1/PD-L1 binding with $IC_{50} = 0.029 \mu\text{g/mL}$. Mouse IgG1, κ was measured as a negative control. (C) Nonbiotinylated cynoPD-L1 competitive binding to cynoPD-1 with $IC_{50} = 7.9 \mu\text{g/mL}$ (152 nM), and Non His-tagged cynoPD-1 competitive binding to cynoPD-L1 with $IC_{50} = 1.22 \mu\text{g/mL}$ (28.6 nM). All IC_{50} values were calculated by using nonlinear regression fitting with GraphPad Prism 7.

Troubleshooting Guide

You will find below recommendations for common situations that you might encounter with your AlphaLISA binding assay. If further assistance is needed, do not hesitate to contact our technical support team for assistance.

Issue	Recommendations and Comments
High background signal	<ul style="list-style-type: none"> • Buffer is not freshly made. Make new. • Incubation time is longer than recommended range.
Low AlphaLISA signal	<ul style="list-style-type: none"> • Optimize EnVision with Plate format.
High variation between replicates or low Z' values	<ul style="list-style-type: none"> • Make sure that reagents are at the bottom of the well by tapping or swirling the plate gently on a smooth surface after each addition.

You will find detailed recommendations for common situations you might encounter with your AlphaLISA Assay kit at:

http://www.perkinelmer.com/in/resources/technicalresources/applicationsupportknowledgebase/alphalisa-alphascreen-no-washassays/alpha_troubleshoot.xhtml

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

This product is not for resale or distribution except by authorized distributors.

LIMITED WARRANTY: PerkinElmer warrants that, at the time of shipment, the above named product is free from defects in material and workmanship and conforms to the specifications set forth above. PerkinElmer makes no other warranty, express or implied with respect to the product and expressly disclaims any warranty of merchantability or fitness for any particular purpose. Notification of any breach of the foregoing warranty must be made within 60 days of receipt of the product, unless otherwise provided in writing by PerkinElmer. No claim shall be honored if the customer fails to notify PerkinElmer within the period specified. The sole and exclusive remedy of the customer for any breach of the foregoing warranty is limited to either the replacement of the non-conforming product or the refund of the invoice price of the product. PERKINELMER SHALL NOT BE LIABLE FOR ANY DIRECT, INDIRECT, SPECIAL, INCIDENTAL, CONSEQUENTIAL OR PUNITIVE DAMAGES, WHETHER BASED ON CONTRACT, TORT, STRICT LIABILITY OR OTHERWISE, ARISING OUT OF THE DESIGN, MANUFACTURE, SALE, DELIVERY, OR USE OF THE PRODUCTS, EVEN IF THE LIMITED REMEDIES PROVIDED HEREIN FAIL OF THEIR ESSENTIAL PURPOSE OR PERKINELMER IS ADVISED OF THE POSSIBILITY OF SUCH DAMAGES.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com

For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.