

Glucagon AlphaLISA Immunoassay Kit

Product number: AL312 HV/C/F

ATTENTION!

Glucagon Analyte included in this kit is provided in solution. Please store at -20°C upon receipt of the kit. Upon the first use of the kit aliquot the analyte and store the aliquots at -20°C to avoid further freeze/thaw cycles.

Caution: For Laboratory Use. A research product for research purposes only.

Lot specific kit information can be found at www.perkinelmer.com/COA

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
General Recommendations.....	5
Assay Procedure.....	5
Data Analysis.....	9
Assay Performance Characteristics.....	9
Troubleshooting Guide.....	11

Product Information

- Application:** This kit is designed for the quantitative determination of Glucagon in cell culture media, sera and plasma, using a homogeneous AlphaLISA assay (no wash steps).
- Sensitivity:** Lower Detection Limit (LDL): 20 pg/mL
Lower Limit of Quantification (LLOQ): 60 pg/mL
EC₅₀: 21 ng/mL
Min/Max counts: 230/ 110 000 counts
- Dynamic range:** 20 - 300 000 pg/mL (Figure 1).

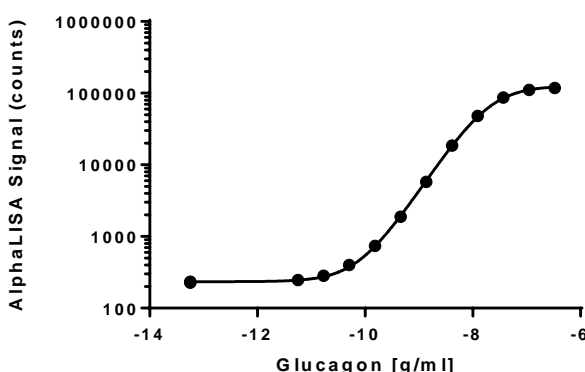


Figure. 1. Typical sensitivity curve in HiBlock buffer. The data was generated using a white Optiplate™-384 microplate and the EnVision® Multilabel Plate Reader with Alpha option 2102.

- Storage:** Immediately store glucagon analyte at -20°C upon arrival upon first use aliquot and store at -20°C to avoid unnecessary freeze/thaw cycles. The remaining kit components must be stored dark at +4°C.
- 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, EC₅₀ and LDL 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 counts may vary between bead lots and the instrument used, with no impact on LDL measurement.

Analyte of Interest

Glucagon is a 29 amino acid peptide hormone produced by the pancreas. Glucagon generally functions as a counter-regulatory hormone opposing the actions of insulin to maintain appropriate levels of blood glucose. Normal human serum glucagon levels range from 50-200 pg/mL. The glucagon:insulin ratio controls the rate of gluconeogenesis and glycogenolysis, disruption of this ratio can have severe metabolic implications. The present kit permits detection of glucagon (i.e. analyte) in different sample matrices.

Description of the AlphaLISA Assay

AlphaLISA technology allows the detection of molecules of interest in buffer, cell culture media, serum and plasma in a highly sensitive, quantitative, reproducible and user-friendly mode. In an AlphaLISA assay, a Biotinylated Anti-Analyte Antibody binds to the Streptavidin-coated Alpha Donor beads, while another Anti-Analyte Antibody is conjugated to AlphaLISA Acceptor beads. In the presence of the analyte, the beads come into close proximity. The excitation of the Donor beads provokes the release of singlet oxygen molecules that triggers a cascade of energy transfer in the Acceptor beads, resulting in a sharp peak of light emission at 615 nm (Figure 2).

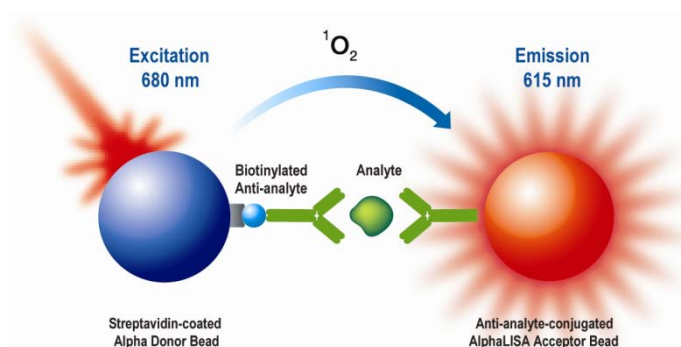


Figure 2. AlphaLISA Assay principle.

Precautions

- The AlphaScreen[®] 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.
- Some analytes are present in saliva. Take precautionary measures to avoid contamination of the reagent solutions.
- The Biotinylated Anti-Analyte Antibody contains sodium azide. Contact with skin or inhalation should be avoided.

Kit Content: Reagents and Materials

Kit components	AL312HV (100 assay points ^{***})	AL312C (500 assay points ^{***})	AL312F (5 000 assay points ^{***})
AlphaLISA Anti-Glucagon Acceptor beads stored in PBS, 0.05% Proclin-300, pH 7.2	20 µL @ 5 mg/mL (1 brown tube, <u>white</u> cap)	50 µL @ 5 mg/mL (1 brown tube, <u>white</u> cap)	500 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% Proclin-300, pH 7.4	100 µL @ 5 mg/mL (1 brown tube, <u>black</u> cap)	200 µL @ 5 mg/mL (1 brown tube, <u>black</u> cap)	200 µL @ 5 mg/mL (2 brown tube, <u>black</u> caps)
Biotinylated Antibody Anti-Glucagon stored in PBS, 0.1% Tween-20, 0.05% NaN ₃ , pH 7.4	100 µL @ 500 nM (1 tube, <u>black</u> cap)	150 µL @ 500 nM (1 tube, <u>black</u> cap)	1.5 mL @ 500 nM (1 tube, <u>black</u> cap)
AlphaLISA Glucagon Analyte Solution (100 µg/mL)	1 tube, <u>clear</u> cap	1 tube, <u>clear</u> cap	1 tube, <u>clear</u> cap
AlphaLISA HiBlock Buffer (10X) **	3 mL, 1 small bottle	10 mL, 1 small bottle	100 mL, 1 large bottle

* The thawed analyte 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. It has been demonstrated that the Glucagon analyte solution is stable for at least 6 months at -20°C. One vial contains an amount of Glucagon sufficient for performing 10 standard curves. Additional vials can be ordered separately (cat # AL312S).

** Contains 250 mM HEPES, pH 7.4, 1% Casein, 10 mg/mL Dextran-500, 5% Triton X-100, 5% BSA and 0.5% Proclin-300. Extra buffer can be ordered separately (cat # AL004C: 10 mL, cat # AL004F: 100 mL).

Note: 10X buffer is slightly tan. If not fully in suspension when diluted to the final 1X solution, it is recommended to centrifuge it for 5 min at 1000 rpm and use the supernatant. It should be noted however, that the appearance of the buffer does not affect its efficacy.

*** The number of assay points is based on an assay volume of 100 µL in 96-well plates (AL312HV) or 50 µL in 96- or 384-well assay plates using the kit components at the recommended concentrations.

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. Note that sodium azide from the Biotinylated Antibody stock solution will not interfere with the AlphaLISA signal (0.0001% final in the assay).

Specific additional required reagents and materials:

The following materials are recommended:

Item	Suggested source	Catalog #
TopSeal™-A Adhesive Sealing Film	PerkinElmer Inc.	6050195
EnVision®-Alpha Reader	PerkinElmer Inc.	-

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 analyte) before use to improve recovery of content (2000g, 10-15 sec). Re-suspend all reagents by vortexing before use.
- Use Milli-Q[®] grade H₂O (18 MΩ•cm) to dilute 10X AlphaLISA HiBlock Buffer to reconstitute the lyophilized analyte.
- When diluting the standard or samples, change tips between each standard or sample dilution. When loading reagents in the assay microplate, change tips between each standard or sample addition and after each set of reagents.
- When reagents are added to the microplate, make sure the liquids are at the bottom of the well.
- Small volumes may be prone to evaporation. It is recommended to cover microplates with TopSeal-A 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.
- The standard curves shown in this technical data sheet are provided for information only. A standard curve must be generated for each experiment. The standard curve should be performed in a similar matrix as the samples (e.g. FBS for serum samples).
- AlphaLISA assays can be performed in cell culture medium with or without phenol red, with the following recommendations: if possible, avoid biotin-containing medium (e.g. RPMI medium) as lower counts and lower sensitivity are expected. Add at least 1% FBS or 0.1% BSA to cell culture medium.

Assay Procedure

IMPORTANT: PLEASE READ THE RECOMMENDATIONS BELOW BEFORE USE

- The protocol described below is an example for generating one standard curve in a 50 µL final assay volume (48 wells, triplicate determinations). The protocols also include testing samples in 452 wells. If a different amount of samples are tested, the volumes of all reagents have to be adjusted accordingly, as shown in the table below. These calculations do not include excess reagent to account for losses during transfer of solutions or dead volumes.
- The standard dilution protocol is provided for information only. As needed, the number of replicates or the range of concentrations covered can be modified.
- When testing plasma or serum samples it is important to add a protease inhibitor immediately upon collection. We recommend 500 KIU of Aprotinin per milliliter of whole blood collected.
- Use of four background points in triplicate (12 wells) is recommended when LDL/LLOQ is calculated. One background point in triplicate (3 wells) can be used when LDL/LLOQ is not calculated.
- Sensitivity can be increased 2 fold (10 pg/ml) by increasing the incubation time of the sample, acceptor beads and biotinylated antibody from 1.5 hours to overnight (12-18 hours). Donor beads should be added after the extended incubation and incubated for 30 minutes at 23° C before reading the assay plate.

		Volume				
Format	# of data points	Final	Sample	AlphaLISA beads / Biotin Antibody MIX	SA-Donor beads	Plate recommendation
AL312HV	100	100 µL	10 µL	40 µL	50 µL	White OptiPlate-96 (cat # 6005290) White ½ AreaPlate-96 (cat # 6005560)
AL312C	250	100 µL	10 µL	40 µL	50 µL	White OptiPlate-96 (cat # 6005290) White ½ AreaPlate-96 (cat # 6005560)
	500	50 µL	5 µL	20 µL	25 µL	White ½ AreaPlate-96 (cat # 6005560) White OptiPlate-384 (cat # 6007290) Light gray AlphaPlate™-384 (cat # 6005350)
	1 250	20 µL	2 µL	8 µL	10 µL	Light gray AlphaPlate-384 (cat # 6005350) ProxiPlate™-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290)
	2 500	10 µL	1 µL	4 µL	5 µL	Light gray AlphaPlate-1536 (cat # 6004350)
AL312F	5 000	50 µL	5 µL	20 µL	25 µL	White ½ AreaPlate-96 (cat # 6005560) White OptiPlate-384 (cat # 6007290) Light gray AlphaPlate-384 (cat # 6005350)
	12 500	20 µL	2 µL	8 µL	10 µL	Light gray AlphaPlate-384 (cat # 6005350) ProxiPlate-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290)
	25 000	10 µL	1 µL	4 µL	5 µL	Light gray AlphaPlate-1536 (cat # 6004350)

High sensitivity protocol (3 incubation steps) – Dilution of standards in 1X AlphaLISA HiBlock Buffer

The protocol described below is recommended when generating one standard curve in a 50 μL final assay volume (48 wells, triplicate determinations with manual pipetting). *If a different amount of samples are tested, the volumes of all reagents have to be adjusted accordingly.*

If higher detection sensitivity is required the sensitivity can be increased 2 fold to 10 $\mu\text{g}/\text{ml}$ by increasing the incubation time of the sample, acceptor beads and biotinylated antibody from 1.5 hours to overnight (12-18 hours). Donor beads should be added after the extended incubation and incubated for 30 minutes at 23° C before reading the assay plate.

Notes: This protocol requires AlphaLISA HiBlock Buffer and it is a 3 step assay!

1) Preparation of 1X AlphaLISA HiBlock Buffer:

- Add 1 mL of 10X AlphaLISA HiBlock Buffer to 9 mL H_2O .

2) Preparation of Glucagon analyte standard dilutions:

- The provided Glucagon solution is provided at 100 $\mu\text{g}/\text{ml}$. The first point of the curve is 1 $\mu\text{g}/\text{ml}$ so a 100 fold dilution of the provided solution is required.
- Prepare standard dilutions as follows (change tip between each standard dilution) in 1X AlphaLISA HiBlock Buffer. **It is not necessary to dilute the standard in the final sample matrix. All spike recovery data provided in this document is interpolated off of a standard curve diluted in the provided HiBlock Buffer.**

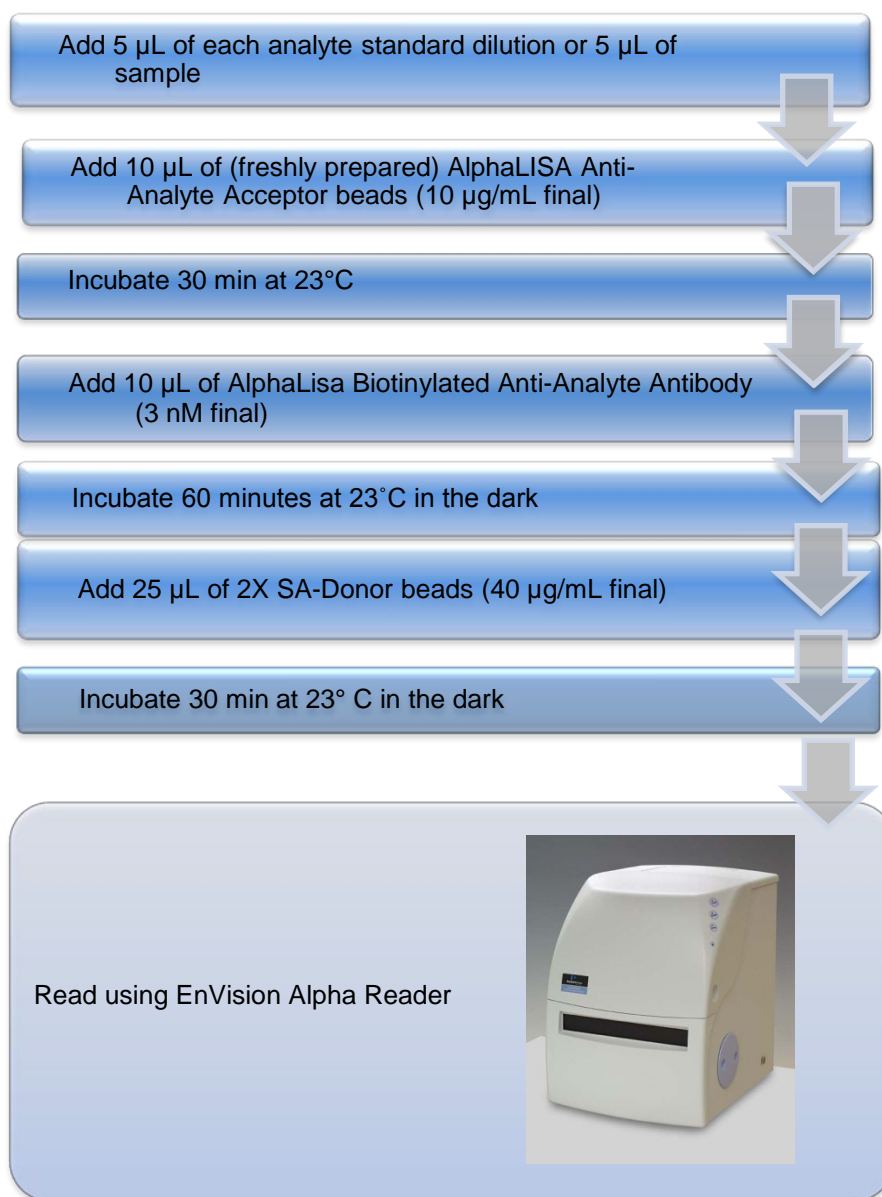
Tube	Vol. of Glucagon (μL)	Vol. of diluent (μL) *	[Glucagon] in standard curve
			($\mu\text{g}/\text{mL}$ in 5 μL)
A	5 μL of reconstituted Glucagon	495	1000000
B	60 μL of tube A	120	333333.3
C	60 μL of tube B	120	111111.1
D	60 μL of tube C	120	37037.0
E	60 μL of tube D	120	12345.7
F	60 μL of tube E	120	4115.2
G	60 μL of tube F	120	1371.7
H	60 μL of tube G	120	457.2
I	60 μL of tube H	120	152.4
J	60 μL of tube I	120	50.8
K	60 μL of tube J	120	16.9
L	60 μL of tube K	120	5.6
M ** (background)	0	100	0
N ** (background)	0	100	0
O ** (background)	0	100	0
P ** (background)	0	100	0

* Dilute standards in diluent (e.g. 1X AlphaLISA HiBlock Buffer).

At low concentrations of analyte, a significant amount of analyte can bind to the vial. Therefore, load the analyte standard dilutions in the assay microplate within 60 minutes of preparation.

** Four background points in triplicate (12 wells) are used when LDL/LLOQ is calculated. If LDL/LLOQ does not need to be calculated, one background point in triplicate can be used (3 wells).

- 3) Preparation of 5X AlphaLISA Anti-Glucagon Acceptor beads (50 µg/mL):
Add 15 µL of 5 mg/mL AlphaLISA Anti-Glucagon Acceptor beads to 1485 µL of 1X AlphaLISA HiBlock Buffer. Prepare just before use.
- 4) Preparation of 5X AlphaLISA Biotinylated Antibody Anti- Glucagon (15 nM):
Add 45 µL of 500 nM Biotinylated Antibody Anti-Glucagon to 1455 µL of 1X AlphaLISA HiBlock Buffer. Prepare just before use.
- 5) Preparation of 2X Streptavidin (SA) Donor beads (80 µg/mL): Keep the beads under subdued laboratory lighting.
Add 48 µL of 5 mg/mL SA-Donor beads to 2952 µL of 1X AlphaLISA HiBlock Buffer.
- 6) Samples:
 - If applicable, dilute samples to be tested in diluent (e.g. 1X AlphaLISA HiBlock Buffer).
- 7) In a 96- or 384-well microplate:



Data Analysis

- Calculate the average count value for the background wells.
- Generate a standard curve by plotting the AlphaLISA counts versus the concentration of analyte. A log scale can be used for either or both axes. No additional data transformation is required.
- Analyze data according to a nonlinear regression using the 4-parameter logistic equation (sigmoidal dose-response curve with variable slope) and a $1/Y^2$ data weighting (the values at maximal concentrations of analyte after the hook point should be removed for correct analysis).
- The LDL is calculated by interpolating the average background counts (12 wells without analyte) + 3 x standard deviation value (average background counts + (3xSD)) on the standard curve.
- The LLOQ as measured here is calculated by interpolating the average background counts (12 wells without analyte) + 10 x standard deviation value (average background counts + (10xSD)) on the standard curve. Alternatively, the true LLOQ can be determined by spiking known concentrations of analyte in the matrix and measuring the percent recovery, and then determining the minimal amount of spiked analyte that can be quantified within a given limit (usually +/- 20% or 30% of the real concentration).
- Read from the standard curve the concentration of analyte contained in the samples.
- If samples have been diluted, the concentration read from the standard curve must be multiplied by the dilution factor.

Assay Performance Characteristics

AlphaLISA assay performance described below was determined using the described protocol.

Sensitivity:

The LDL and LLOQ were calculated as described above. The values correspond to the lowest concentration of analyte that can be detected in a volume of 5 μ L using the recommended assay conditions.

LDL (pg/mL)	LLOQ (pg/ml)	Buffer/Media used	# of experiments
20	60	AlphaLISA HiBlock Buffer	8

* Note that LDL/ LLOQ can be decreased (i.e. sensitivity increased) by increasing the volume of analyte in the assay (e.g. use 10 μ L of analyte in a final assay volume of 50 μ L).

Assay precision:

The following assay precision data were calculated from the eight independent assays using two different kit lots. In each lot, the analytes were prepared in AlphaLISA HiBlock Buffer (HBB). Each assay consisted of one standard curve comprising 12 data points (each in triplicate) and 12 background wells (no analytes). The assays were performed in 384-well format using AlphaLISA HiBlock Buffer.

- Intra-assay precision:

The intra-assay precision was determined using a total of 6 independent determinations in triplicate. Shown CV%.

Glucagon (ng/ml)	HBB
1.4	12%

- Inter-assay precision:

The inter-assay precision was determined using a total of 6 independent determinations.

Glucagon (ng/ml)	HBB
1.4	13%

- Recovery:

Three known concentrations of analyte were spiked into cell culture media containing 10% FBS, cell culture media containing 1% BSA, depleted human sera and AlphaLISA HiBlock Buffer (HBB). All samples were run alongside a standard curve diluted in AlphaLISA HiBlock Buffer, this standard curve was used to interpolate the concentrations of the samples. The percent recovery is defined as assay measured concentration with respect to the spiked concentration. The average recovery from two independent measurements is reported.

Spike (Glucagon ng/mL)	% Recovery			
	AlphaLISA HiBlock Buffer	DMEM with FBS	DMEM with BSA	Depleted Human Sera
10	115	124	124	110
15	92	121	127	117
1	87	103	101	115

Troubleshooting Guide

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 BioSignal Inc. warrants that, at the time of shipment, the products sold by it are free from defects in material and workmanship and conform to specifications which accompany the product. PerkinElmer BioSignal Inc. makes no other warranty, express or implied with respect to the products, including any warranty of merchantability or fitness for any particular purpose. Notification of any breach of warranty must be made within 60 days of receipt unless otherwise provided in writing by PerkinElmer BioSignal Inc. No claim shall be honored if the customer fails to notify PerkinElmer BioSignal Inc. within the period specified. The sole and exclusive remedy of the customer for any liability of PerkinElmer BioSignal Inc. of any kind including liability based upon warranty (express or implied whether contained herein or elsewhere), strict liability contract or otherwise is limited to the replacement of the goods or the refunds of the invoice price of goods. PerkinElmer BioSignal Inc. shall not in any case be liable for special, incidental or consequential damages of any kind.

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

