

## HTRF LCL161 Red Ligand

**Part # 64BDIAPB3RED**

**Test Size#:** 10 Binding assays (K<sub>d</sub> determination) - assay volume: 20 µL

**Revision:** #03 of September 2023      **Store at:** ≤-16°C

**This product is intended for research purposes only. It is not intended to be used for therapeutic or diagnostic purposes.**

### ASSAY PRINCIPLE

Revvity' HTRF LCL161 Red Ligand (#Revvity 64BDIAPB3RED) is primarily intended to perform affinity binding curves using HTRF® technology and can be used in association with the HTRF cIAP1 BIR3 Binding kit (#Revvity 64BDIAPB3PEG/H).

The HTRF LCL161 Red Ligand is particularly suited for cooperativity binding studies to assess the effect of a PROTAC protein substrate on the affinity of the LCL161 Red Ligand for cIAP1 BIR3 protein. HTRF LCL161 Red Ligand binding is detected in a direct binding assay format using an anti-GST antibody labeled with Terbium Cryptate which binds to Human cIAP1 BIR3 GST-tagged. When the dyes are in close proximity, the excitation of the donor with a light source (laser or flash lamp) triggers a Fluorescence Resonance Energy Transfer (FRET) towards the acceptor, which in turn fluoresces at a specific wavelength (665 nm). The specific binding signal is calculated by subtracting the non-specific binding signal from the total signal, enabling determination of K<sub>d</sub> for HTRF LCL161 Red Ligand (Fig. 1)

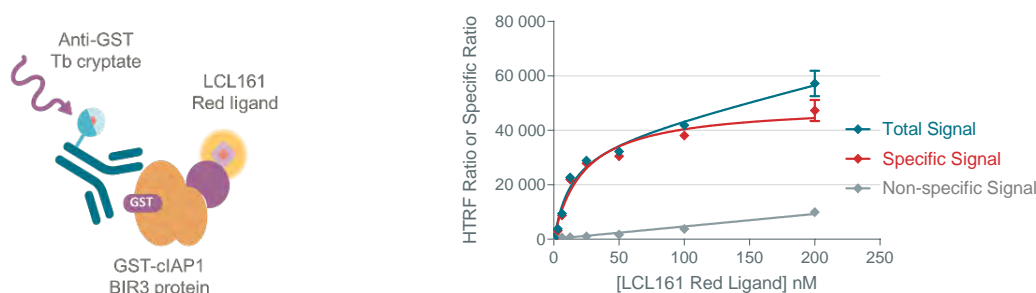
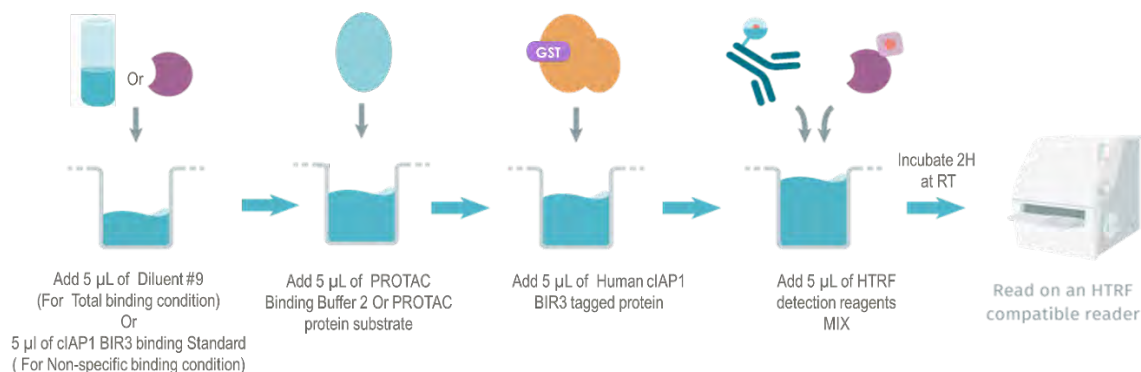


Figure 1: Principle of K<sub>d</sub> determination with HTRF LCL161 Red Ligand.

### MANUAL AT A GLANCE



Assay volumes can be adjusted proportionally to run the assay in 96 or 1536 well microplates.

Make sure you use the appropriate setup for Terbium Cryptate. For more information about setup and HTRF® compatible readers, please visit our website at: [www.revvity.com](http://www.revvity.com)

**MATERIAL PROVIDED:**

HTRF LCL161 Red Ligand - 5X	1 vial - 200 µL Frozen	
-----------------------------	---------------------------	---

**PURCHASE SEPARATELY:**

- cIAP1 BIR3 Standard (#Revvity 64BDIAPB3CDA)
- HTRF PROTAC Binding Buffer 2 (#Revvity 64BDE32RDF)
- Diluent 9 (5X) (#Revvity 62DL9DDA/C)
- HTRF cIAP1 BIR3 Binding kit (#Revvity 64BDIAPB3PEG/H)
- Low volume white (only) microplate\*
- HTRF®-Certified Reader \*\*. Make sure the setup for Terbium Cryptate is used.

\* For HTRF microplate recommendations, please visit [www.revvity.com](http://www.revvity.com)

\*\* For a list of HTRF-compatible readers and setup recommendations, please visit [www.revvity.com](http://www.revvity.com)

**STORAGE AND STABILITY**

Store the HTRF LCL161 Red Ligand at -16°C or below. Under appropriate storage conditions, HTRF LCL161 Red Ligand is stable until the expiry date indicated on the label.

To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -16°C or below. Aliquot and store rapidly.

Volume of HTRF LCL161 Red Ligand aliquots should not be under 10 µL.

Once thawed, aliquoted HTRF LCL161 Red Ligand can be frozen once.

Working solutions should be prepared the day of the experiment and not stored for longer period.

**REAGENT PREPARATION****BEFORE YOU BEGIN:**

- It is recommended to prepare all reagents in the specified buffers. The use of incorrect buffers may affect reagent stability and assay results.
- cIAP1 BIR3 binding standard will be used to determine the non-specific binding signal.
- PROTAC Binding Buffer 2 is ready-to-use.
- cIAP1 BIR3 GST tagged protein must be prepared following instructions from the package insert of the HTRF cIAP1 BIR3 Binding Kit.
- Diluent 9 (5X) must be homogenized with a vortex then diluted 5-fold with distilled water (e.g. 200µL of Diluent 9 (5X) + 800 µL of distilled water) and mixed gently to obtain the Diluent 9 (1X).

**► cIAP1 BIR3 BINDING STANDARD**

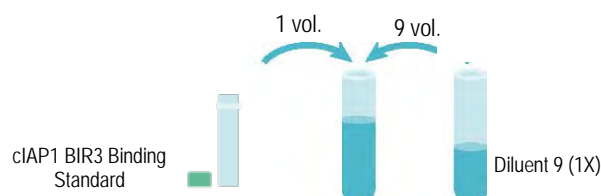
- To prepare the cIAP1 BIR3 binding standard Stock solution:

Thaw the cIAP1 BIR3 binding Standard. Centrifuge.

The stock solution (100 µM) can be frozen and stored at -16°C or below.

- To prepare the cIAP1 BIR3 binding standard working solution:

Homogenize the cIAP1 BIR3 binding Standard stock solution then dilute it 10-fold with Diluent 9 (1X) (e.g. 20µL of cIAP1 BIR3 binding Standard stock solution + 180µL Diluent 9 (1X)). Mix gently. cIAP1 BIR3 binding standard working concentration is 10 µM.



### ► LCL161 RED LIGAND

- To prepare the LCL161 Red Ligand Stock solution:

Thaw the HTRF LCL161-Red Ligand. Centrifuge.

This 5X stock solution can be frozen and stored at -16°C or below.

- To prepare the LCL161 Red Ligand working solutions:

Procedure to prepare one binding curve of the HTRF LCL161-Red Ligand (for monitoring Total and Non-specific Binding):

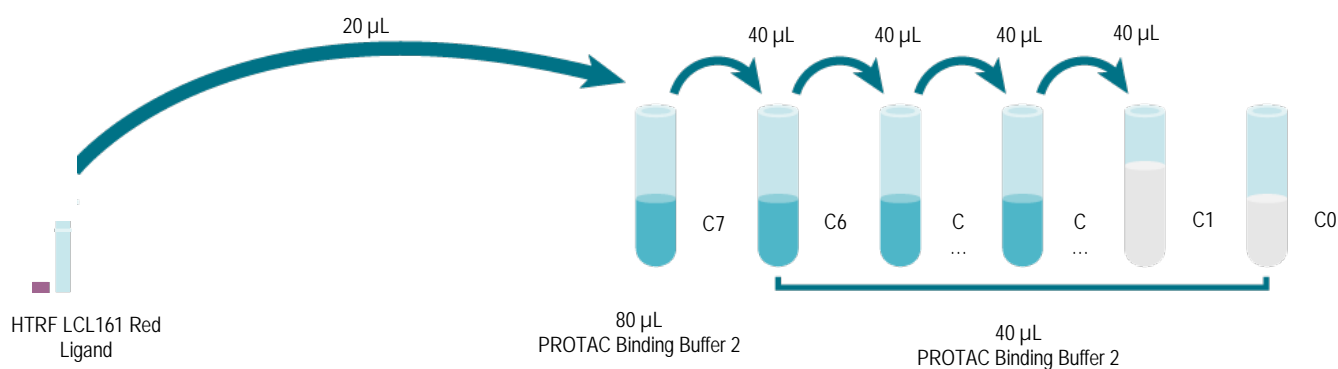
Dilute the LCL161 Red Ligand stock solution 5-fold with PROTAC Binding Buffer 2 to prepare high LCL161-Red Ligand (C7), (e.g. 20  $\mu$ L of LCL161-Red Ligand stock solution + 80  $\mu$ L of PROTAC Binding Buffer 2). Mix gently.

Use the high LCL161-Red Ligand (C7) to prepare LCL161-Red Ligand recommended range of concentrations using 1/2 serial dilutions as follows:

- Dispense 40  $\mu$ L of PROTAC Binding Buffer 2 into each vial from C6 to C1.
- Add 40  $\mu$ L of LCL161 Red Ligand to 40  $\mu$ L of PROTAC Binding Buffer 2, mix gently and repeat the 1/2 serial dilution to make standard solutions: C6, C5, C4, C3, C2, C1.
- C0 is PROTAC Binding Buffer 2 buffer alone (Negative control).

This will create 7 concentrations for the LCL161 Red Ligand to be used for Total and Non-specific Binding conditions.

*In order to counteract any sticking, we recommend changing tips between each dilution.*



### ► GST TB CRYPTATE ANTIBODY:

- To prepare the GST Tb Cryptate Stock solution:

Thaw the GST Tb Cryptate antibody included in the cIAP1 BIR3 Binding kit (#Revvity 64BDIAPB3PEG/H). Centrifuge.

This stock solution can be frozen and stored at -16°C or below.

- To prepare the GST Tb Cryptate working solution:

Homogenize the GST Tb Cryptate antibody stock solution then dilute it 25-fold with PROTAC Binding Buffer 2, (e.g. 80  $\mu$ L of GST Tb Cryptate antibody stock solution + 1920  $\mu$ L of PROTAC Binding Buffer 2). Mix gently.

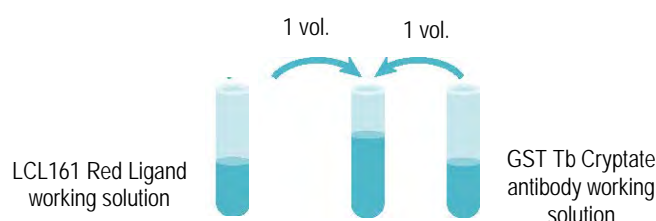


**WARNING:** Do not prepare the mAb anti-GST Tb Cryptate working solution following instructions from the package insert of the cIPA1 BIR3 Binding kit (#Revvity 64BDIAPB3PEG/H).

► **DETECTION REAGENT MIX SOLUTIONS OF HTRF LCL161-RED LIGAND + GST TB CRYPTATE ANTIBODY:**







Just prior dispensing, mix each working solution of HTRF LCL161 Red Ligand and GST Tb Cryptate antibody by adding 1 volume of GST Tb Cryptate antibody working solution to 1 volume of each HTRF LCL161 Red Ligand working solutions (C7 to C0) (e.g. 50  $\mu$ L of HTRF LCL161 Red Ligand working solution + 50  $\mu$ L of GST Tb Cryptate antibody working solution).

This will create 7 Mix for the LCL161 Red Ligand and GST Tb Cryptate antibody.



MIX	MIX PREPARATION	LCL161 RED LIGAND CONCENTRATION IN THE MIX SOLUTION (nM)	LCL161 RED LIGAND FINAL CONCENTRATION (nM)
Mix 7	50 $\mu$ L C7 LCL161-Red Ligand + 50 $\mu$ L of working solution of GST Tb Cryptate antibody	800	200
Mix 6	50 $\mu$ L C6 LCL161-Red Ligand + 50 $\mu$ L of working solution of GST Tb Cryptate antibody	400	100
Mix 5	50 $\mu$ L C5 LCL161-Red Ligand + 50 $\mu$ L of working solution of GST Tb Cryptate antibody	200	50
Mix 4	50 $\mu$ L C4 LCL161-Red Ligand + 50 $\mu$ L of working solution of GST Tb Cryptate antibody	100	25
Mix 3	50 $\mu$ L C3 LCL161-Red Ligand + 50 $\mu$ L of working solution of GST Tb Cryptate antibody	50	12.5
Mix 2	50 $\mu$ L C2 LCL161-Red Ligand + 50 $\mu$ L of working solution of GST Tb Cryptate antibody	25	6.25
Mix 1	50 $\mu$ L C1 LCL161-Red Ligand + 50 $\mu$ L of working solution of GST Tb Cryptate antibody	12.5	3.125
Mix 0	50 $\mu$ L PROTAC Binding Buffer 2 + 50 $\mu$ L of working solution of GST Tb Cryptate antibody	0	0

## ASSAY MANUAL

	NON-SPECIFIC BINDING SIGNAL	TOTAL BINDING SIGNAL
<b>Step 1</b> 	Dispense 5 $\mu$ L of cIAP1 BIR3 Binding Standard into each well	Dispense 5 $\mu$ L of Diluent#9 (1X) into each well
<b>Step 2</b> 	Add 5 $\mu$ L of PROTAC Binding Buffer 2 or 5 $\mu$ L of PROTAC substrate protein (diluted in the PROTAC Binding Buffer 2)	
<b>Step 3</b> 	Add 5 $\mu$ L of Human cIAP1 BIR3 GST Tagged protein	
<b>Step 4</b> 	Add 5 $\mu$ L of premixed HTRF LCL161 Red Ligand and GST Tb Cryptate antibody working solutions to all wells	
<b>Step 5</b> 	Seal the plate and incubate 2 hours at RT	
<b>Step 6</b> 	Remove the plate sealer and read on an HTRF® compatible reader	

## EXAMPLE OF MAP PLATE

	NON-SPECIFIC BINDING			TOTAL BINDING								
	1	2	3	4				5	6			
A	5 µL of cIAP1 BIR3 Binding Standard	Repeat Well A1	Repeat Well A1	5 µL of Diluent #9 (1X) 5 µL of PROTAC Binding Buffer 2 or PROTAC protein substrate 5 µl Human cIAP1 BIR3 GST Tagged protein 5 µL MIX 0 (PROTAC Binding Buffer 2 and GST Tb Cryptate antibody)				Repeat Well A4	Repeat Well A4			
B	5 µL of cIAP1 BIR3 Binding Standard	Repeat Well B1	Repeat Well B1	5 µL of Diluent #9 (1X) 5 µL of PROTAC Binding Buffer 2 or PROTAC protein substrate 5 µl Human cIAP1 BIR3 GST Tagged protein 5 µL MIX 1 (LCL161-Red Ligand C1 and GST Tb Cryptate antibody)				Repeat Well B4	Repeat Well B4			
C	5 µL of cIAP1 BIR3 Binding Standard	Repeat Well C1	Repeat Well C1	5 µL of Diluent #9 (1X) 5 µL of PROTAC Binding Buffer 2 or PROTAC protein substrate 5 µl Human cIAP1 BIR3 GST Tagged protein 5 µL MIX 2 (LCL161-Red Ligand C2 and GST Tb Cryptate antibody)				Repeat Well C4	Repeat Well C4			
D	5 µL of cIAP1 BIR3 Binding Standard	Repeat Well D1	Repeat Well D1	5 µL of Diluent #9 (1X) 5 µL of PROTAC Binding Buffer 2 or PROTAC protein substrate 5 µl Human cIAP1 BIR3 GST Tagged protein 5 µL MIX 3 (LCL161-Red Ligand C3 and GST Tb Cryptate antibody)				Repeat Well D4	Repeat Well D4			
E	5 µL of cIAP1 BIR3 Binding Standard	Repeat Well E1	Repeat Well E1	5 µL of Diluent #9 (1X) 5 µL of PROTAC Binding Buffer 2 or PROTAC protein substrate 5 µl Human cIAP1 BIR3 GST Tagged protein 5 µL MIX 4 (LCL161-Red Ligand C4 and GST Tb Cryptate antibody)				Repeat Well E4	Repeat Well E4			
F	5 µL of cIAP1 BIR3 Binding Standard	Repeat Well F1	Repeat Well F1	5 µL of Diluent #9 (1X) 5 µL of PROTAC Binding Buffer 2 or PROTAC protein substrate 5 µl Human cIAP1 BIR3 GST Tagged protein 5 µL MIX 5 (LCL161-Red Ligand C5 and GST Tb Cryptate antibody)				Repeat Well F4	Repeat Well F4			
G	5 µL of cIAP1 BIR3 Binding Standard	Repeat Well G1	Repeat Well G1	5 µL of Diluent #9 (1X) 5 µL of PROTAC Binding Buffer 2 or PROTAC protein substrate 5 µl Human cIAP1 BIR3 GST Tagged protein 5 µL MIX 6 (LCL161-Red Ligand C6 and GST Tb Cryptate antibody)				Repeat Well G4	Repeat Well G4			
H	5 µL of cIAP1 BIR3 Binding Standard	Repeat Well H1	Repeat Well H1	5 µL of Diluent #9 (1X) 5 µL of PROTAC Binding Buffer 2 or PROTAC protein substrate 5 µl Human cIAP1 BIR3 GST Tagged protein 5 µL MIX 7 (LCL161-Red Ligand C7 and GST Tb Cryptate antibody)				Repeat Well H4	Repeat Well H4			

A	1	2	3	4	5	6	7	8	9	10	11	12

	1	2	3	4	5	6	7	8	9	10	11	12
A												
B												
C												
D												
E												
F												
G												
H												
I												

## DATA REDUCTION & INTERPRETATION

1. Calculate the ratio of the acceptor and donor emission signals for each individual well.

$$\text{Ratio} = \frac{\text{Signal 665 nm}}{\text{Signal 620 nm}} \times 10^4$$

2. Calculate the % CVs. The mean and standard deviation can then be worked out from ratio replicates.

$$\text{CV (\%)} = \frac{\text{Standard deviation}}{\text{Mean Ratio}} \times 100$$

3. Calculate the specific Binding signal for each concentration used of HTRF LCL161 Red Ligand.

$$\text{Specific Binding signal} = \text{Total Binding signal} - \text{Non-specific Binding signal}$$

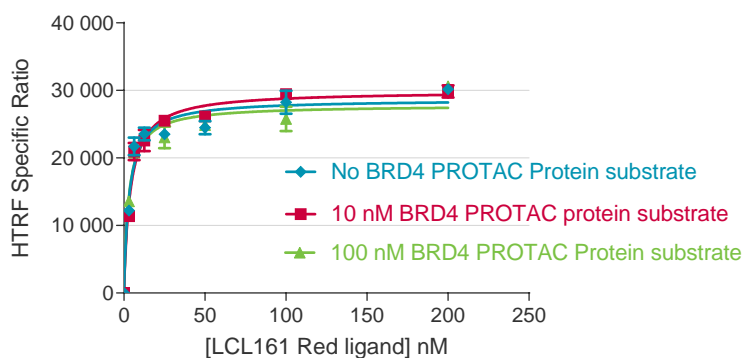
For more information about data reduction, please visit [www.revvity.com](http://www.revvity.com)

## RESULTS

This data must not be substituted for the data obtained in the laboratory and should be considered only as an example (readouts on an HTRF compatible reader). Results may vary from one HTRF® compatible reader to another.

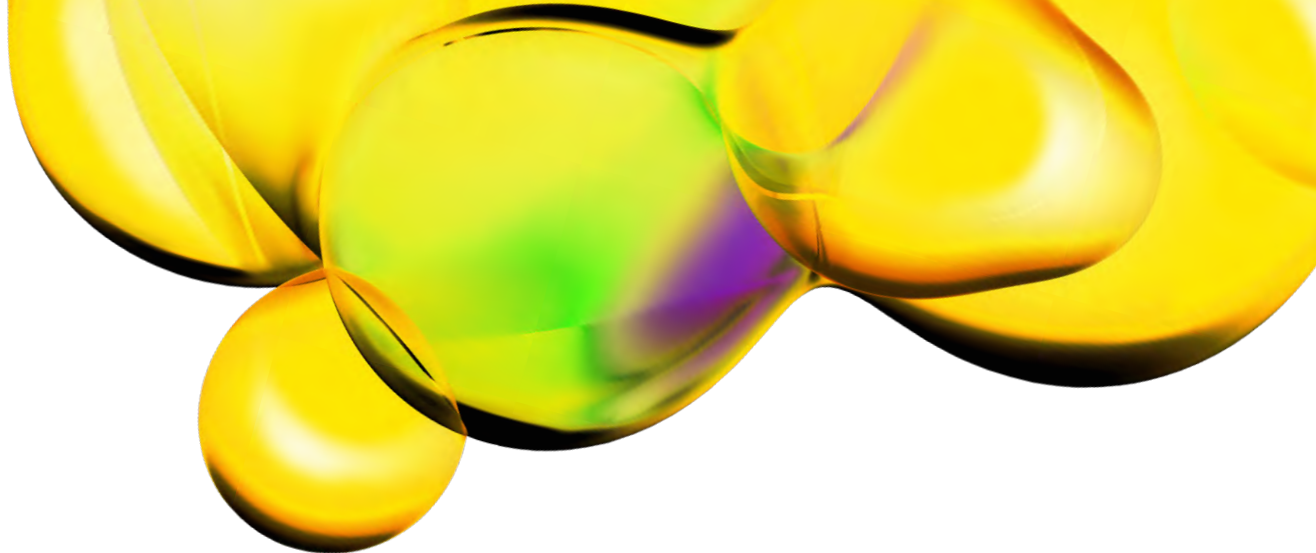
In this example, the affinity binding (K<sub>d</sub>) of the LCL161-Red Ligand to GST-cIAP1 BIR3 protein was assessed in the absence or presence of the targeted protein BRD4 (10nM or 100nM). This result indicates that the presence of BRD4 does not change the K<sub>d</sub> of the Red Ligand.

Consequently, this example shows that a cooperativity experiment with a PROTAC molecule composed of a cIAP ligand and a BRD4 warhead can be set up. Finally, the alpha factor can be established by dividing the K<sub>i</sub> of the PROTAC in the binary complex (cIAPBIR3 -PROTAC) by the K<sub>i</sub> of the PROTAC in the ternary complex (cIAPBIR3 -PROTAC-BRD4).



## ANALYTICAL CHARACTERISTICS

HTRF LCL 161 Red Ligand K <sub>d</sub> (reference without PROTAC protein substrate)	12 nM ±10 (2SD)
---	-----------------



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

Manufactured by Cisbio Bioassays - Parc Marcel Boiteux - 30200 Codolet - FRANCE

[www.revvity.com](http://www.revvity.com)

revvity

**Revvity, Inc.**  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.