



## PolyGT-Biotin

Part # 61GT0BLA and 61GT0BLD

Test size: 5,000 tests (61GT0BLA), 100,000 tests (61GT0BLD) - assay volume: 20  $\mu$ L

Revision: #10 of September 2023

Store at 2-8°C

For research use only. Not for use in therapeutic or diagnostic procedures.

### REAGENT DESCRIPTION

Poly-(Glu,Tyr) (4:1) is a random copolymer which has been shown to be a substrate for tyrosine-specific protein kinases. Poly-(GT)-biotin may be used with Streptavidin-XL665 and Anti-phosphotyrosine-cryptate for easily optimized HTRF<sup>®</sup> tyrosine kinase assays.

### MATERIALS

Reagent	5,000 tests	100,000 tests
PolyGT-Biotin Lyophilized - MW: 44kDa	1 vial	1 vial

Revvity reagents - Not provided	Part #
HTRF KinEASE detection buffer 200 mL - ready-to-use	62SDBRDF
Plates - HTRF 96-well low volume plate	66PL96001

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

For reading, an HTRF<sup>®</sup>-certified reader is needed. Make sure to use the setup for Eu<sup>3+</sup> Cryptate.

For a list of HTRF<sup>®</sup>-certified readers and setup recommendations, please visit [www.revvity.com](http://www.revvity.com)

### STORAGE AND STABILITY

- Store the lyophilized reagent at 2-8°C.
- Under appropriate storage conditions, reagents are stable until the expiry date indicated on the batch information.
- Upon reconstitution, the reagent stock solution is stable 2 weeks at 2-8°C. It can be refrozen (at  $\leq$ -16°C) and thawed at least two more times.

### ASSAY FORMAT

When used as suggested, one vial will provide sufficient reagent for 5,000 tests and 100,000 tests respectively on the basis of 50 ng of Poly-(GT) per well (1.14 pmol/well) using a 20  $\mu$ L final assay volume.

## REAGENT HANDLING

### BUFFERS

Revvity KinEASE detection buffer (#62SDBRDF) has been optimized for maximum performance and is ready to use.

When using specific in-house buffers for the preparation of working solutions, make sure to use a **phosphate-free buffer** (i.e. hepes), with a **pH maintained between 5.5 and 8.5**. It can be supplemented with BSA (0.1%), and detergents such as Tween 20, Triton X100 and CHAPS (up to 0.5%) to prevent reagent coating. Avoid SDS, due to its denaturing effect on XL665.

Use of Europium antibody conjugate solution requires a final KF concentration between 100 mM and 400 mM.

It is mandatory to use the same buffer to prepare the donor and the acceptor (d2 or XL665) conjugates.

### CONJUGATES

Allow the lyophilized reagent to warm up at room temperature for at least 30 minutes.

Prepare the stock & working solutions according to the table below

PolyGT-Biotin*	Stock solution preparation	Working solution preparation
5,000 tests	Reconstitute with 250 $\mu$ L distilled water. Mix gently.	Dilute the stock solution in KinEASE detection buffer. Mix gently.
100,000 tests	Reconstitute with 5 mL distilled water. Mix gently.	Dilute the stock solution in KinEASE detection buffer. Mix gently.

\*Additional info is included on the batch information provided with the reagent

After reconstitution, the stock solution can be divided into aliquots and frozen for additional use.

Make sure to prepare stock and working solutions according to the instructions that correspond to the packaging you have purchased (number of tests).

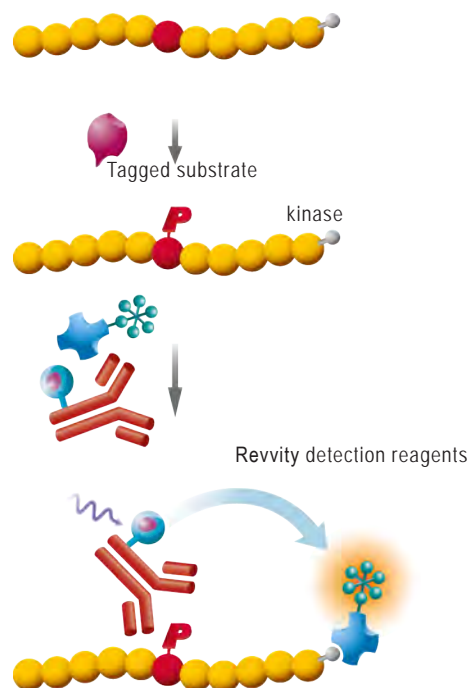
## COMPANION REAGENTS

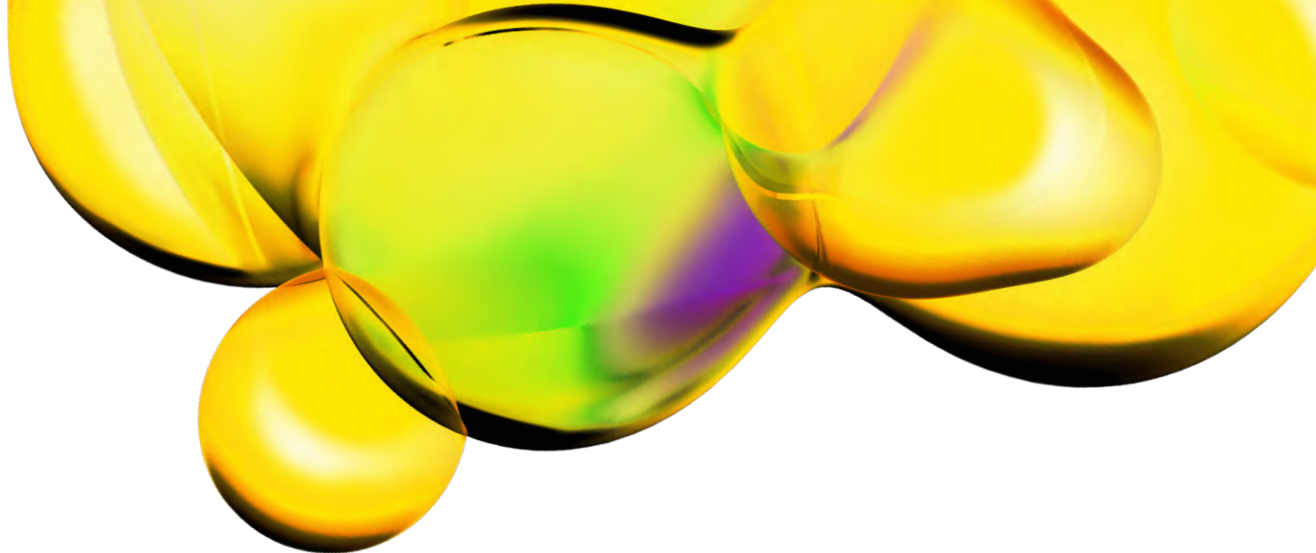
As illustrated beside, all kinase assays are based on the same format.

The enzymatic reaction is usually carried out with a biotinylated substrate (protein, peptide and the enzyme itself in the case of autophosphorylation).

The phosphorylated substrate is then detected using the specific anti phosphoresidue antibody coupled to  $\text{Eu}^{3+}$  Cryptate and a XL665 conjugate such as streptavidin-XL665<sup>entl</sup> (ref 611SAXLA) or Streptavidin-XL665 (ref 610SAXLA).

Alternatively, other tags such as GST-, 6HIS, c-myc or DNP may be used instead of biotin to label the kinase substrate.





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.