



PAb Anti Human IgG-Tb cryptate

Part # 61HFCTAF, 61HFCTAA & 61HFCTAB

Test size: 1,000 tests (61HFCTAF), 5,000 tests (61HFCTAA), 20,000 tests (61HFCTAB)

Assay volume: 20 µL

Revision: #05 of September 2023

Store at ≤ -16°C

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

REAGENT DESCRIPTION

In an HTRF protein:protein interaction assay, one protein is labeled (directly or indirectly) with the donor, and the other protein is labeled (directly or indirectly) with the acceptor. When the two proteins interact, the donor molecule is brought into proximity with the acceptor molecule. Excitation of the donor will result in signal generation proportional to the binding of proteins.

IgGs from goat anti-Human immunoserum were isolated by immunoaffinity chromatography using Human IgG coupled to agarose beads. The purified polyclonal antibodies have been labeled with Lumi4® Tb. This reagent may be used to detect the binding of unlabeled human specific antibody on a target molecule or on Ig fusion proteins.

This reagent may cross-react with immunoglobulins from other species. Therefore, it is always recommended to control the appropriate blanks levels in all types of assays specially when IgG from other animal species are involved

MATERIALS

Reagent	1,000 tests	5,000 tests	20,000 tests
PAb Anti Human IgG-Tb cryptate Frozen in phosphate buffer pH 7.0 containing protease free bovine serum albumin	1 vial - 0.05 mL	1 vial - 0,25 mL	1 vial - 1 mL

Revvity reagents Not provided	Part #
PPI - Terbium detection buffer	Cat # 61DB10RDF* *200 mL - ready-to-use
Plates - HTRF 96-well low volume plate	66PL96001

For HTRF microplate recommendations, please visit www.revvity.com

For reading, an HTRF® - compatible reader is needed. Make sure to use the appropriate set-up.

For a list of HTRF® - compatible readers and set-up recommendations, please visit www.revvity.com

STORAGE AND STABILITY

- Store the reagent at ≤ -16°C.
- Under appropriate storage conditions, reagents are stable until the expiry date indicated on the label.
- Once thawed, stock solutions are stable for two days at 2-8°C. They can be refrozen (at ≤ -16°C) and thawed once only. Do not repeat freezing and thawing.

ASSAY FORMAT

When used as suggested, one vial from the two available sizes will provide sufficient reagent for 1,000 tests, 5,000 tests and 20,000 tests respectively, using a 20 μ L final assay volume.

	Volume
Other assay components	10 μ L
Acceptor (d2 or XL665) conjugate	5 μ L
Donor (Eu or Tb Cryptate) conjugate	5 μ L
Final volume*	20 μ L

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

REAGENT HANDLING

BUFFERS

Revvity PPI - Terbium detection buffer Cat # 61DB10RDF* has been optimized for maximum performance.

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

When using specific in-house buffers for the preparation of working solutions, we recommend a basic buffer such as PBS (Phosphate Buffered Saline) or 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 or CHAPS (up to 0.5%) to prevent reagent coating. Avoid SDS, due to its denaturing effect on XL665.

Please note that a phosphate-free buffer must be used in biochemical kinase assays to prevent interferences pertaining to the binding of anti-phospho specific antibodies.

Use of Terbium conjugates does not require KF.

CONJUGATES

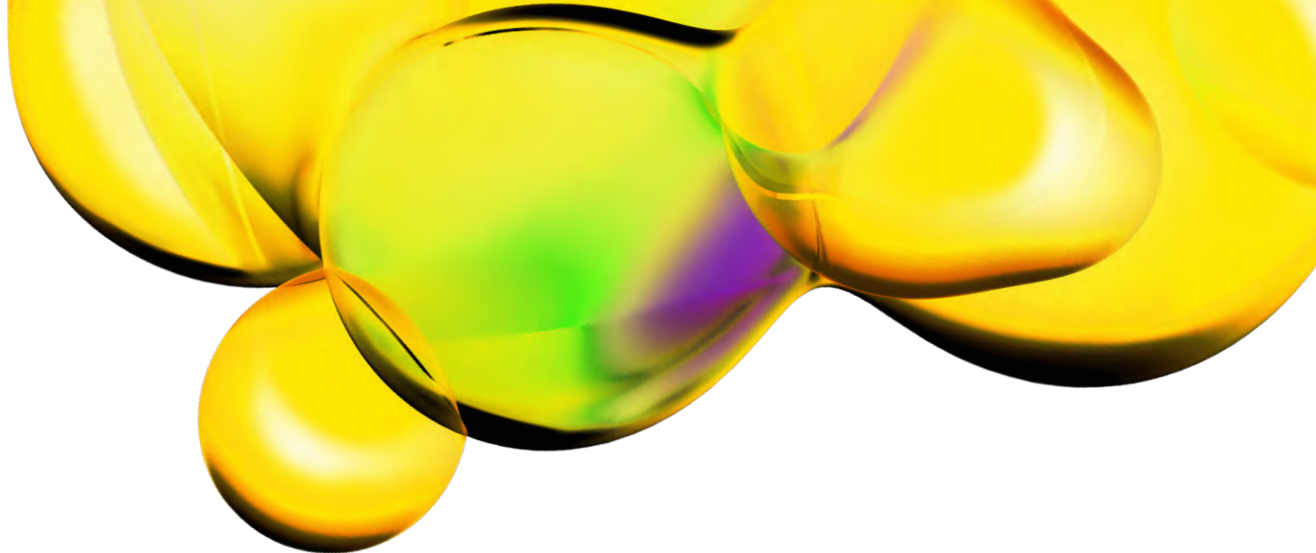
Allow each vial of frozen conjugate to warm up at room temperature.

Mix gently.

PAb Anti Human IgG-Tb cryptate	Stock solution preparation	Working solution preparation (see assay format above)
1,000 tests	0.05 mL Mix gently	Dilute 100-fold the stock solution in PPI detection buffer Mix gently E.g. Add 4.95 mL PPI detection buffer to 0.05 mL of stock solution
5,000 tests	0.25 mL Mix gently	Dilute 100-fold the stock solution in PPI detection buffer Mix gently E.g. Add 24.75 mL PPI detection buffer to 0.25 mL of stock solution
20,000 tests	1 mL Mix gently	Dilute 100-fold the stock solution in PPI detection buffer Mix gently E.g. Add 99 mL PPI detection buffer to 1 mL of stock solution

Additional info is included in the batch information provided with the reagent.

The optimal amount per well will be dependent on assay conditions. For additional information on assay optimization, please refer to the technical note: Guidelines for optimizing protein : protein interaction assays using HTRF PPI reagents. Make sure to prepare stock and working solutions according to the instructions that correspond to the packaging you have purchased (number of tests).



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

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.