



MAb PY20-Eu cryptate

PROTOCOL

Part # 61Y20KLA and 61Y20KLB

Test size: 5,000 tests (61Y20KLA), 20,000 tests (61Y20KLB) - assay volume: 20 μ L

Revision: 09 (Dec. 2017)

Store at 2-8°C

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

REAGENT DESCRIPTION

PY20 is a mouse monoclonal IgG2b which was produced using Phosphotyrosine conjugated to BSA as immunogen. This monoclonal was shown to react specifically with phosphorylated tyrosine, both as free amino acid or when conjugated to carriers. No cross-reactivity has been observed with nonphosphorylated tyrosine, phosphothreonine, phosphoserine, AMP or ATP. It is suitable for use in a wide variety of tyrosine kinase activity tests involving receptor or non-receptor enzymes, and specific or ubiquitous substrates.

MATERIALS

Reagent	5,000 tests	20,000 tests
MAb PY20-Eu cryptate Lyophilized	1 vial	1 vial

Cisbio 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 <http://www.cisbio.com/microplate-recommendations>.

For reading, an HTRF[®]-certified reader is needed. Make sure to use the setup for Eu³⁺ Cryptate.

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

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 1 week at 2-8°C. It can be refrozen (at \leq -16°C) and thawed once only.

ASSAY FORMAT

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

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

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

REAGENT HANDLING

BUFFERS

Cisbio 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

MAB PY20-Eu cryptate*	Stock solution preparation	Working solution preparation (see assay format)
5,000 tests	Reconstitute with 250 μ L distilled water. Mix gently.	Dilute 100-fold the stock solution in KinEASE detection buffer. Mix gently. E.g. Add 24.75 mL of KinEASE detection buffer to 0.25 mL of stock solution.
20,000 tests	Reconstitute with 1 mL distilled water. Mix gently.	Dilute 100-fold the stock solution in KinEASE detection buffer. Mix gently. E.g. Add 99 mL of KinEASE detection buffer to 1 mL of stock solution.

*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).

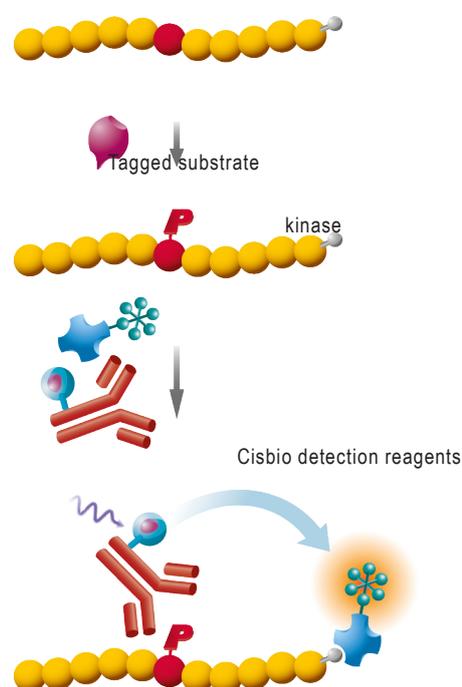
COMPANION REAGENTS

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 Eu^{3+} Cryptate and a XL665 conjugate such as streptavidin-XL^{ent} (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.



This product contains material of biological origin. Use for research purposes only. Do not use in humans or for diagnostic purposes. The purchaser assumes all risk and responsibility concerning reception, handling and storage.

Copyright 2017 Cisbio, France. All rights reserved. HTRF, Tag-lite, and EPIgeneous are trademarks or registered trademarks of Cisbio. All other trademarks are the property of their respective owners.

FOR MORE INFORMATION

Europe and other countries +33(0)466-796-705 U.S. and Canada 1-888-963-4567 China +86-21-5018-9880
Japan +81-(0)43-306-8712 Visit www.cisbio.com to find a list of our regional distributors