



STREPTAVIDIN-EU CRYPTATE KINASE BINDING

Part # 62KBSAKAF & 62KBSAKAB

Test size: 1,000 tests (# 62KBSAKAF), 20,000 tests (# 62KBSAKAB)

Assay volume: 20µL

Form: Frozen

Store at: -16°C or below

Revision: #02 of September 2023

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

REAGENT DESCRIPTION

The Streptavidin-Eu cryptate Kinase Binding is used in the HTRF® Kinase Binding format in combination with a N-terminal Biotinylated Kinase and red-fluorescent tracers.

MATERIALS

Reagent	1,000 tests	20,000 tests
Streptavidin-Eu cryptate Kinase Binding 100 X	1 vial - 0.05 mL	1 vial - 1 mL

Revvity reagents - NOT PROVIDED	Amount & Part #
Kinase Binding Buffer	20 mL - Part # 62KBBRDD 200 mL - Part# 62KBBRDF
Staurosporine-Red	1 nmol - Part# 62KB01REDC 20 nmol - Part# 62KB01REDE
Dasatinib-Red	1 nmol - Part# 62KB02REDC 20 nmol - Part# 62KB02REDE
Sunitinib-Red	1 nmol - Part# 62KB03REDC 20 nmol - Part# 62KB03REDE

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

** For a list of HTRF-compatible readers and setup recommendations, please visit www.revvity.com

STORAGE AND STABILITY

- Store the reagent at ≤ -16°C.
- Once thawed, the solution can be frozen once (≤ -16°C or below).
- Do not repeat freeze/thaw cycles. To avoid them, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -16°C or below.

ASSAY FORMAT

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

For additional information, please refer to the [HTRF® Kinase Binding Guide](#).

Saturation Kinase Binding Assay K_D determination		Competition Kinase Binding Assay IC_{50} - K_I determination	
Reagent	Volume	Reagent	Volume
Kinase Binding Buffer	5 μ L	Inhibitor	5 μ L
Tagged-kinase	5 μ L	Tagged-kinase	5 μ L
Streptavidin-Eu cryptate Kinase Binding	5 μ L	Streptavidin-Eu cryptate Kinase Binding	5 μ L
Inhibitor-Red	5 μ L	Inhibitor-Red	5 μ L
Final volume	20 μ L	Final volume	20 μ L

REAGENT HANDLING

BUFFER

The Revvity Kinase Binding Buffer (Part# 62KBBRDD & Part#/62KBBRDF) has been optimized for maximum performance.

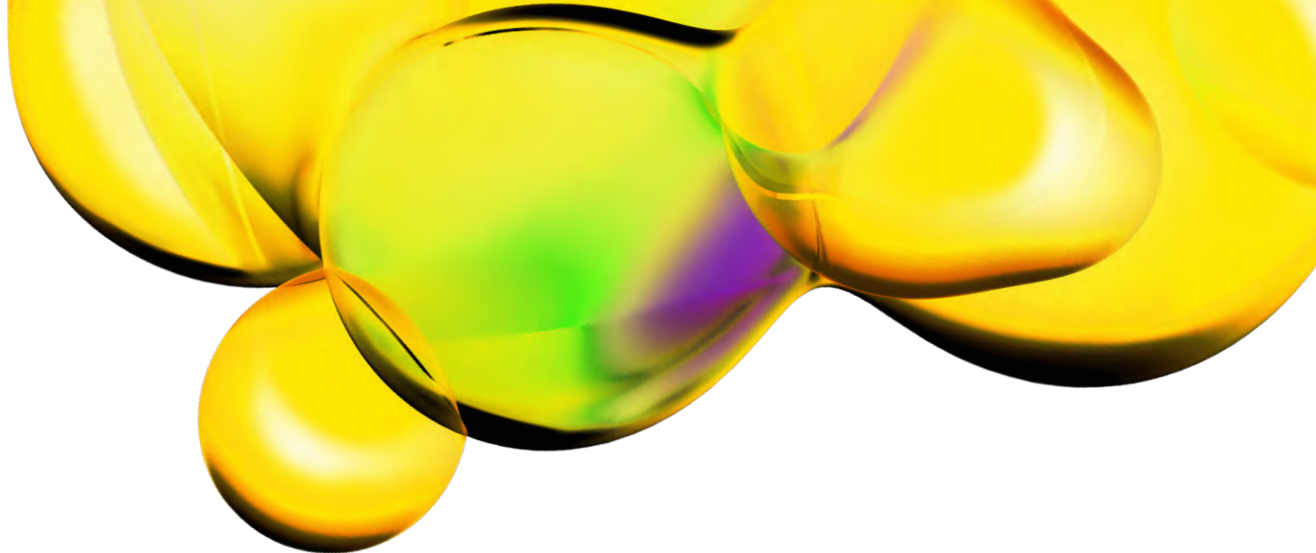
The use of an incorrect diluent may affect reagent stability and assay results.

We recommend filtering (0.22 μ m) the buffer before use.

STREPTAVIDIN-EU CRYPTATE KINASE BINDING

Before use, allow the reagent to warm up at room temperature and homogenize it with a vortex.

Streptavidin-Eu cryptate Kinase Binding 100X	Working solution preparation
1,000 tests - 0.05 mL	Dilute 100-fold the stock solution in Kinase Binding Buffer Mix gently. e.g. 0.05 mL of stock solution + 4.95 mL Kinase Binding Buffer
20,000 tests - 1 mL	Dilute 100-fold the stock solution in Kinase Binding Buffer Mix gently. e.g. 1 mL of stock solution + 99 mL Kinase Binding Buffer



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.