



MAB ANTI GST-EU CRYPTATE KINASE BINDING

PROTOCOL

Part # 62KBGSTKAF & 62KBGSTKAB

Test size: 1,000 tests (# 62KBGSTKAF), 20,000 tests (# 62KBGSTKAB)
Assay volume: 20 μ L

Form: Frozen

Store at: -16°C or below

Revision: 01 - March 2019

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



REAGENT DESCRIPTION

The MAb Anti GST-Eu cryptate Kinase Binding is used in the HTRF® Kinase Binding format in combination with a GST-tagged Kinase and red-fluorescent tracers.

MATERIALS

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

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

** For a list of HTRF-compatible readers and setup recommendations, please visit <http://www.cisbio.com/compatible-readers>

STORAGE AND STABILITY

- Store the reagent at \leq -16°C.
- Once thawed, the solution can be frozen once (\leq -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
MAB Anti GST-Eu cryptate Kinase Binding	5 μ L	MAB Anti GST-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 Cisbio 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.

MAB ANTI GST-EU CRYPTATE KINASE BINDING

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

MAB Anti GST-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

This product contains material of biologic 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 2019 Cisbio. All rights reserved.

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