

HTRF® Research Reagents

HTRF Kinase-6HIS Bind. Discovery kit

Product Number : 62KBD02PEA Lot Number : 01K

Expiration date : 2025-02-19 Document version : 1

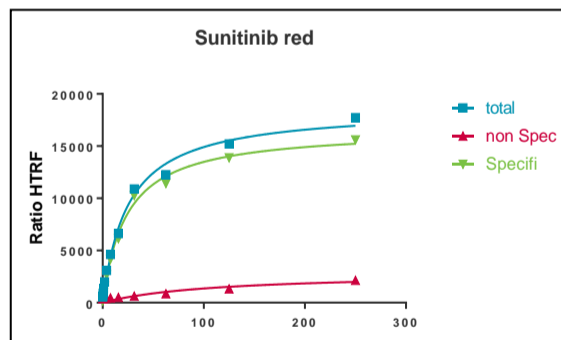
Reagents :	Lot n°
HTRF Dasatinib-Red - 0.5 nmol	04RA
HTRF Kinase Bind. Buf. - 20 mL	04RA
HTRF Staurosporine-Red - 0.5 nmol	04RA
HTRF Sunitinib-Red - 0.5 nmol	04RA
HTRF Anti 6HIS-Eu - Kinase Bind. 1K pts	01RA

Results :

Incubation : 1h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : One site binding (hyperbola)

Assay characteristics :

	Found value	Specifications Rev 01
KD	25	20-50 nM
S/B	8,3	> 6



HTRF Kinase-6HIS Bind. Discovery kit - Lot n° : 01K

Accepted

Rejected

Quality control date : 2025-02-21	Supervisor's initials : MP
-----------------------------------	----------------------------

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. 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.