

HTRF® Research Reagents

HTRF KRAS G12C/SOS1 PPI kit - 10K pts

Product Number : 64KRASG12PEH Lot Number : 04B

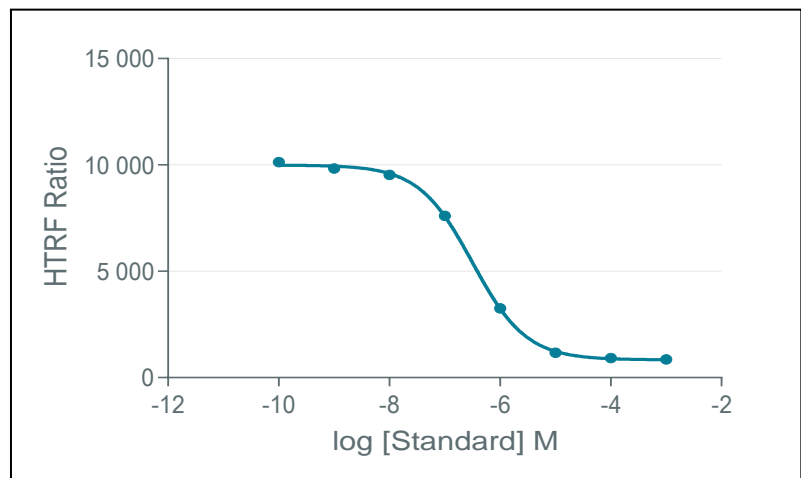
Expiration date : 2025-02-24 Document version : 1

Reagents :	Lot n°
HTRF KRAS G12C/SOS1 PPI kit - 10K pts Tag2 Human SOS1	04RA
HTRF KRAS G12C/SOS1 PPI kit - 10K pts Anti Tag2 Tb Cryptate antibody	03RA
HTRF KRAS G12C/SOS1 PPI kit - 10K pts Tag1 Human KRAS G12C	04RA
HTRF KRAS G12C/SOS1 PPI kit - 10K pts Anti-tag1 XL665 antibody	04RA
HTRF KRAS WT/SOS1 PPI kit - 10K pts GTP	03RA
HTRF KRAS/SOS1 PPI kit - Std	04RA
HTRF Bind. Domain - Detect. Buf.2-130 mL	14RA

Results :

Incubation : 2h at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : Pherastar Plus  
 Fitting model : log(inhibitor) vs.response-Variable slope ( four parameters)

STD	Conc. (µMf)	RATIO
Std 0	0	10137
Std 1	0,001	9837
Std 2	0,01	9530
Std 3	0,1	7611
Std 4	1	3255
Std 5	10	1174
Std 6	100	924
Std 7	1000	861



Assay characteristics :

	Found value	Specifications Rev 01
S/B	14	12-16
IC 50	318	180-700 nM

HTRF KRAS G12C/SOS1 PPI kit - 10K pts - Lot n° : 04B

Accepted

Rejected

Quality control date : 2024-01-26	Supervisor's initials : FL
-----------------------------------	----------------------------

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.