

HTRF® Research Reagents

HTRF KRAS WT/SOS1 PPI kit - 500 pts

Product Number : 64KRASWTPEG Lot Number : 03A

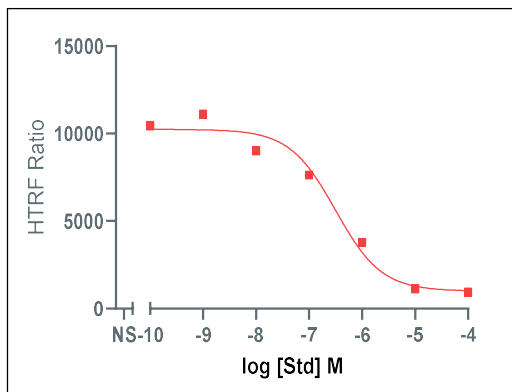
Expiration date : 2025-10-05 Document version : 1

Reagents :	Lot n°
HTRF KRAS WT/SOS1 PPI kit - 500 pts Tag2 Human SOS1	05A
HTRF KRAS/SOS1 PPI kit - Std	04AR
HTRF KRAS WT/SOS1 PPI kit - 500 pts GTP	03RA
HTRF KRAS WT/SOS1 PPI kit - 500 pts Anti Tag2 Tb Cryptate antibody	03A
HTRF KRAS WT/SOS1 PPI kit - 500 pts Tag1 Human KRAS WT	05A
HTRF KRAS WT/SOS1 PPI kit - 500 pts Anti-tag1 XL665 antibody	02RA
HTRF Bind. Domain - Detect. Buf. 2- 20 mL	11RA

Results :

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

STD	Conc. (µMf)	RATIO
Std 0	0	10 445
Std 1	0,001	11 119
Std 2	0,01	9 000
Std 3	0,1	7 627
Std 4	1	3 779
Std 5	10	1 137
Std 6	100	944
Std 7	1000	922



Assay characteristics :

	Found value	Specifications Rev 01
S/B	13	11-15
IC 50	315	300-600 nM

HTRF KRAS WT/SOS1 PPI kit - 500 pts - Lot n° : 02E

Accepted

Rejected

Quality control date : 2024-01-16	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.