

HTRF® Research Reagents

HTRF KRAS WT/SOS1 PPI kit - 500 pts

Product Number : 64KRASWTPEG Lot Number : 03C

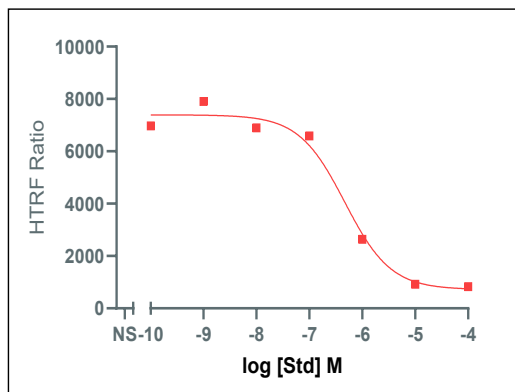
Expiration date : 2027-02-24 Document version : 1

Reagents :		Lot n°
HTRF KRAS WT/SOS1 PPI kit - 500 pts Tag2 Human SOS1		05A
HTRF KRAS/SOS1 PPI kit - Std		05A
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		11B
/		/

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	6 989
Std 1	0,001	7 906
Std 2	0,01	6 900
Std 3	0,1	6 591
Std 4	1	2 635
Std 5	10	917
Std 6	100	829
Std 7	1000	814



Assay characteristics :

	Found value	Specifications Rev 01
S/B	11	11-15
IC 50	504	300-600 nM

HTRF KRAS WT/SOS1 PPI kit - 500 pts - Lot n° : 03C

☒ Accepted☐ Rejected

Quality control date :	2025-08-06	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.