

HTRF® Research Reagents

HTRF KRAS G12C/SOS1 PPI kit - 500 pts

Product Number : 64KRASG12PEG

Lot Number : 03K

Expiration date : 2026-06-27

Document version :

Reagents :		Lot n°
HTRF KRAS G12C/SOS1 PPI kit - 500 pts Tag2 Human SOS1		05A
HTRF KRAS G12C/SOS1 PPI kit - 500 pts Anti Tag2 Tb Cryptate antibody		03RA
HTRF KRAS G12C/SOS1 PPI kit - 500 pts Tag1 Human KRAS G12C		05A
HTRF KRAS G12C/SOS1 PPI kit - 500 pts Anti-tag1 XL665 antibody		03RA
HTRF KRAS/SOS1 PPI kit - Std		05A
HTRF KRAS WT/SOS1 PPI kit - 500 pts GTP		03RA
HTRF Bind. Domain - Detect. Buf. 2- 20 mL		11B
/		/

Results :

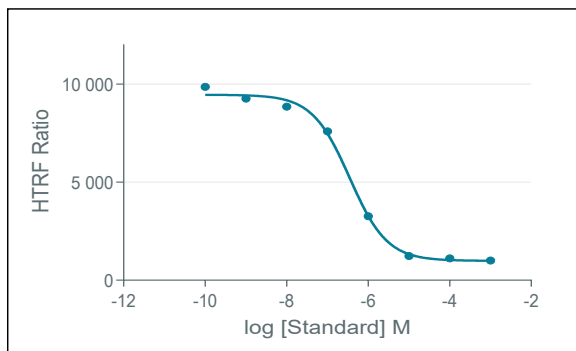
Incubation : 2h at RT

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	9866
Std 1	0,001	9268
Std 2	0,01	8850
Std 3	0,1	7594
Std 4	1	3261
Std 5	10	1232
Std 6	100	1118
Std 7	1000	1008



Assay characteristics :

	Found value	Specifications Rev 01
S/B	13	12-16
IC 50	351	180-700 nM

HTRF KRAS G12C/SOS1 PPI kit - 500 pts - Lot n° : 03K

☒ Accepted☐ Rejected

Quality control date : 2025-08-11

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.