

HTRF® Research Reagents

HTRF KRAS G12C GTP Bind. kit - 500 pts

Product Number : 64BDKRASGPEG Lot Number : 02A

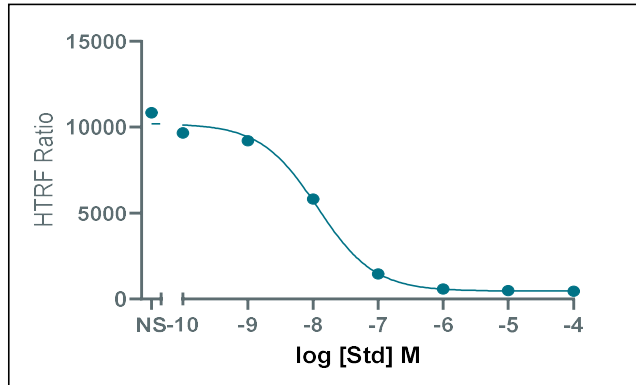
Expiration date : 2025-11-28 Document version : 1

Reagents :	Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL	07RA
HTRF KRAS GTP Bind. kit - Std	01RA
HTRF KRAS G12C GTP Bind. kit - 500 pts Eu Cryptate antibody	01RA
HTRF KRAS G12C GTP Bind. kit - 500 pts Human KRAS G12C 6His-tagged protein	03A
HTRF KRAS G12C GTP Bind. kit - 500 pts GTP-Red reagent	02A

Results :

Incubation : 1h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : /

STD	Conc. (µMf)	RATIO
Std 0	0	10838
Std 1	0,0001	9668
Std 2	0,001	9203
Std 3	0,01	5816
Std 4	0,1	1450
Std 5	1	576
Std 6	10	489
Std 7	100	461



Assay characteristics :

	Found value	Specifications Rev 01
S/B	24	17-24
IC 50	12	12 - 25 nM

HTRF KRAS G12C GTP Bind. kit - 500 pts - Lot n° : 02A

Accepted

Rejected

Quality control date : 2024-08-06	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.