

HTRF® Research Reagents

HTRF (h) RIG-I Bind. Kit - 500 pts

Product Number : 64BDRIGIPEG Lot Number : 04D

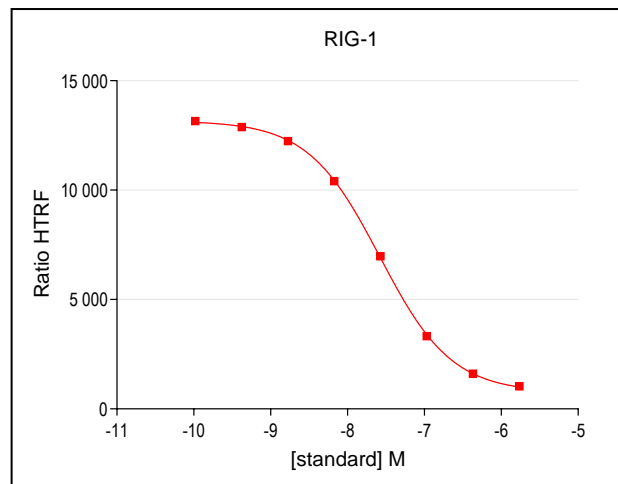
Expiration date : 2025-12-18 Document version : 1

Reagents :	Lot n°
HTRF Diluent 13 - 8 mL	01RA
HTRF Detect. Buf. 16 - 2 mL	01RB
HTRF (h) RIG-I Bind. Kit - Std	04RA
HTRF (h) RIG-I Bind. Kit - 500 pts Streptavidin d2	04AR
HTRF (h) RIG-I Bind. Kit - 500 pts Biot-dsRNA	02RB
HTRF (h) RIG-I Bind. Kit - 500 pts Tb Cryptate antibody	04AR
HTRF (h) RIG-I Bind. Kit - 500 pts protein RIG-His	03RB

Results :

Incubation : 1h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

3P -RNA	Conc. (nMi)	RATIO
Std 0	0	13158
Std 1	0,42	12878
Std 2	1,7	12239
Std 3	6,7	10408
Std 4	27	6979
Std 5	108	3322
Std 6	430	1609
Std 7	1720	1029



Assay characteristics :

	Found value	Specifications Rev 01
IC 50	26	13 - 36 nM
S/B	13	10 - 13

HTRF (h) RIG-I Bind. Kit - 500 pts - Lot n° : 04D

Accepted

Rejected

Quality control date : 2024-02-02	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.