

HTRF® Research Reagents

HTRF TOM20 Kit - 500 pts

Product Number : 64TOM20PEG Lot Number : 01D

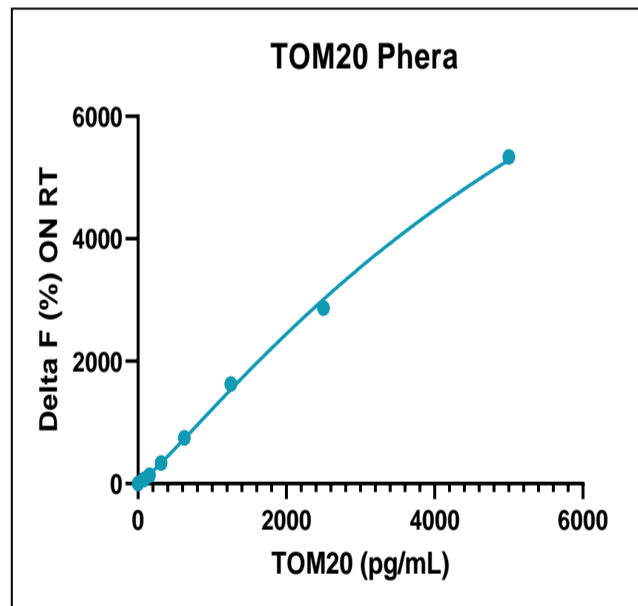
Expiration date : 2024-06-22 Document version : 1

Reagents :	Lot n°
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF TOM20 Kit - Std	01RA
HTRF TOM20 Kit - 500 pts d2 antibody	01RA
HTRF TOM20 Kit - 500 pts Eu Cryptate antibody	01RA
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	12RJ
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	26RA

Results :

Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : 4 Parameter Logistic (4PL)

TOM 20	Conc (ng/ml)	Delta F %	Delta F% Accepted Range
Std 0	0	0%	0-0
Std 1	78	65%	28-78
Std 2	156	139%	117-142
Std 3	312	339%	280-348
Std 4	625	748%	688-765
Std 5	1250	1626%	1549-1622
Std 6	2500	2869%	2637-3385
Std 7	5000	5335%	4813-5471



HTRF TOM20 Kit - 500 pts - Lot n° : 01D

Accepted Rejected

Quality control date : 2023-12-21	Supervisor's initials : FMV
-----------------------------------	-----------------------------

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.