

HTRF® Research Reagents

HTRF Human IL18 kit - 500 pts

Product Number : 62HIL18PEG Lot Number : 06E

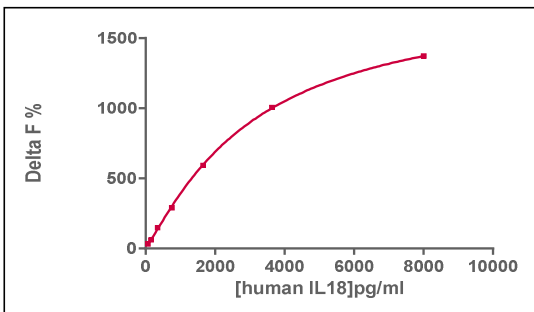
Expiration date : 2025-01-22 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	19G
HTRF Detect. Buf. 3 - 1.5 mL	28A
HTRF Human IL18 kit - Std	03RA
HTRF Human IL18 kit - 500 pts d2 antibody	06RA
HTRF Human IL18 kit - 500 pts Eu Cryptate antibody	06RA
HTRF Human IL18 kit - modop	7

Results :

Incubation : 1h at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : Pherastar Plus  
 Fitting model : 4 Parameter Logistic (4PL 1/y2)\* model

hIL18	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	70.6	33%	22-35
Std 2	155	60%	47-70
Std 3	341	148%	104-155
Std 4	751	291%	228-321
Std 5	1 653	593%	446-652
Std 6	3 636	1006%	845-1097
Std 7	8 000	1372%	1227-1592



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	8.6 pg/ml	≤ 70.6 pg/ml

HTRF Human IL18 kit - 500 pts - Lot n° : 06E

Accepted

Rejected

Quality control date :	2023-11-29	Supervisor's initials :	PL
------------------------	------------	-------------------------	----

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.