

HTRF® Research Reagents

# HTRF Human IL8 kit - 500 pts

Product Number : 62HIL08PEG Lot Number : 06D

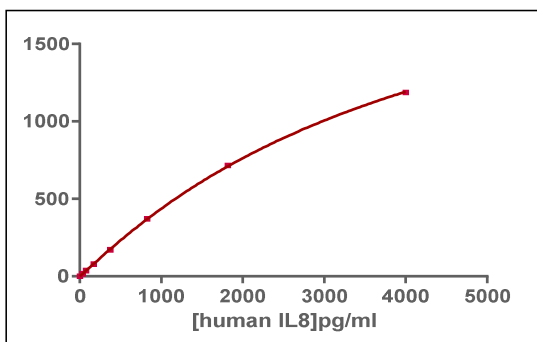
Expiration date : 2025-10-05 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	28RA
HTRF Human IL8 kit - Std	04RA
HTRF Human IL8 kit - 500 pts d2 antibody	06RA
HTRF Human IL8 kit - 500 pts Eu Cryptate antibody	06RA
HTRF Human IL8 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

hIL8	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	35.3	15%	11-22
Std 2	77.6	37%	23-46
Std 3	170.75	79%	55-99
Std 4	375.7	171%	119-213
Std 5	826.4	370%	241-476
Std 6	1818.2	716%	501-978
Std 7	4000	1187%	938-1862



Assay characteristics :

	Found value	Specifications Rev 02
Detection limite	7 pg/ml	≤ 35.3 pg/ml

HTRF Human IL8 kit - 500 pts - Lot n° : 06E

Accepted

Rejected

Quality control date : 2024-01-03	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.