

HTRF® Research Reagents

HTRF Human GM-CSF kit - 500 pts

Product Number :62HGMCSFPEG

Lot Number :06B

Expiration date : 2027-01-23

Document version :1

Reagents :		Lot n°
HTRF Diluent 5 (5X) -2 mL		21B
HTRF Detect. Buf. 3 - 1,5 mL		30B
HTRF Human GM-CSF kit - 500 pts d2 antibody		06A
HTRF Human GM-CSF kit - 500 pts Eu Cryptate antibody		06A
HTRF Human GM-CSF kit - Std		06A

Results :

Incubation :3h at RT

Plate :384 wells small volume white plate 20 µL

Reader :Pherastar Plus

Fitting model :4 Parameter Logistic (4PL 1/γ2)* model

hGM-CSF	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	75	31%	21-38
Std 2	164,9	72%	57-76
Std 3	362,9	148%	125-160
Std 4	798,3	306%	277-322
Std 5	1756,2	559%	539-609
Std 6	3863,6	957%	948-972
Std 7	8500	1277%	1161-1227

* This batch has been validated despite Std 7 out of range

Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	10 pg/ml	≤ 75 pg/ml

HTRF Human GM-CSF kit - 500 pts - Lot n° : 06B

☒ Accepted

☐ Rejected

Quality control date :2024-07-01	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.