

HTRF® Research Reagents

HTRF Human IL5 kit - 500 pts

Product Number : 62HIL05PEG Lot Number : 05A

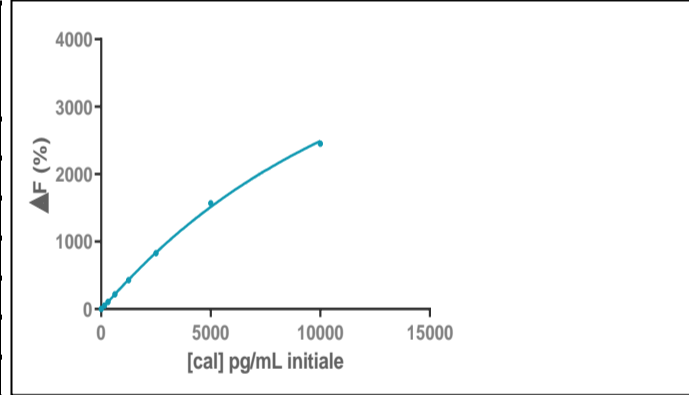
Expiration date : 2025-04-18 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF Human IL5 kit - Std	03RA
HTRF Human IL5 kit - 500 pts Eu Cryptate antibody	05A
HTRF Human IL5 kit - 500 pts XL665 antibody	05A

Results :

Incubation : 2h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : 4 Parameter Logistic (4PL 1/y2)* model

hIL5	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	156	53%	44-84
Std 2	312	106%	84-157
Std 3	625	218%	171-296
Std 4	1250	429%	343-537
Std 5	2500	827%	619-944
Std 6	5000	1568%	1144-1646
Std 7	10000	2401%	1747-2410



Assay characteristics :

	Found value	Specifications Rev 01
Detection limite	11,8 pg/ml	≤ 156 pg/ml

HTRF Human IL5 kit - 500 pts - Lot n° : 05A

Accepted

Rejected

Quality control date : 2024-04-24	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.