

HTRF® Research Reagents

HTRF CHO HCP Kit - 500 pts

Product Number : 64CHOPEG Lot Number : 02A

Expiration date : 2025-06-12 Document version : 1

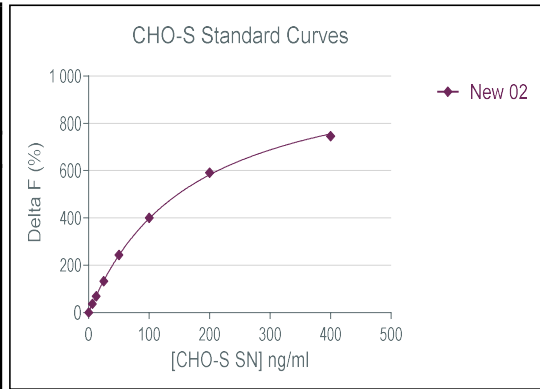
Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 17 - 1.5 mL	01RA
HTRF CHO HCP Kit - Std	01RB
HTRF CHO HCP Kit - 500 pts d2 antibody	02A
HTRF CHO HCP Kit - 500 pts Eu Cryptate antibody	02A

Results :

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

Assay characteristics :

CHO	Conc (ng/ml)	Delta F %	Delta F% Accepted Range Provisoire
Std 0	0	0%	0-0
Std 1	6.25	37%	25 - 37
Std 2	12.5	69%	51 - 77
Std 3	25	133%	99 - 149
Std 4	50	244%	182 - 273
Std 5	100	400%	314 - 470
Std 6	200	591%	484 - 727
Std 7	400	745%	630 - 946



HTRF CHO HCP Kit - 500 pts - Lot n° : 02A

Accepted

Rejected

Quality control date : 2024-02-09	Supervisor's initials : AC
-----------------------------------	----------------------------

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.