

HTRF® Research Reagents

HTRF Mouse IL6 kit - 500 pts

Product Number : 62MIL06PEG Lot Number : 08B

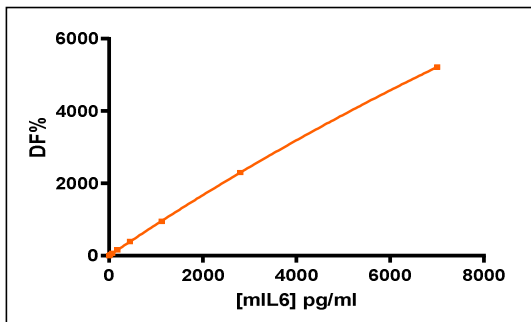
Expiration date : 2025-05-06 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF Mouse IL6 kit - Std	04RA
HTRF Mouse IL6 kit - 500 pts d2 antibody	08RA
HTRF Mouse IL6 kit - 500 pts Eu Cryptate antibody	08RA
HTRF Mouse IL6 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

M IL6	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	28.7	33%	24-43
Std 2	71.7	72%	49-112
Std 3	179.2	162%	137-252
Std 4	448	392%	373-619
Std 5	1120	953%	891-1471
Std 6	2800	2300%	2160-3281
Std 7	7000	5217%	4899-5814



Assay characteristics :

	Found value	Specifications Rev 01
Detection limite	3.5 pg/ml	≤ 28.7 pg/ml

HTRF Mouse IL6 kit - 500 pts - Lot n° : 08B

Accepted

Rejected

Quality control date : 2023-11-17	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.