

HTRF® Research Reagents

HTRF Human SMN kit - 500 pts

Product Number : 63ADK033PEG Lot Number : 14D

Expiration date : 2026-11-29 Document version : 1

Reagents :	Lot n°
HTRF Detect. Buf. 3 - 7 mL	30AR
HTRF Human SMN kit - 500 pts d2 antibody	14B
HTRF Human SMN kit - 500 pts Eu Cryptate antibody	14B
HTRF Human SMN kit - Ctrl lysate	14A
HTRF P-T prot. - Lysis Buf.2 (4X) 2 mL	13AR

Results :

 Incubation : Over night at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Envision
 Fitting model : /

Assay characteristics :

	Found value	Specifications Rev 01
Negative	409	
Control Lysate	1924	
S/B	4,5	> 4

HTRF Human SMN kit - 500 pts - Lot n° : 14D

 Accepted

 Rejected

Quality control date : 2024-05-24	Supervisor's initials : FL
-----------------------------------	----------------------------

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.