

HTRF® Research Reagents

Biotin-Heparan sulfate-Eu cryptate Control

Product number : 61BHSKAA Lot Number : 25RB

Manufacturing date : 25/07/2018 Document version : 1

EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing enzyme concentrations (Heparanase).

Biotin-Heparan sulfate-Eu is frozen in PO4 buffer 100 mM pH7 ; 0.1% Tween.

ASSAY FORMAT

Biotin-Heparan sulfate-Eu-Cryptate + Streptavidin-d2 + Heparanase

REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
Streptavidin-d2	19A	2.5 µg/mL
Biotin-Heparan sulfate-Eu-Cryptate	24A	202.5ng/mL
Heparanase	INSIGHT 08/2013	12.5-> 1 000 ng/mL

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Biotin-Heparan sulfate-Eu-Cryptate current batch	25RB	202.5ng/mL

Other reagents were the same as those used for the reference test.

EXPERIMENTAL PROCEDURE

Reagents were thawed, diluted and dispensed as follows:

- 5 µL Biotin-HS-Cryptate
- 5 µL Heparanase

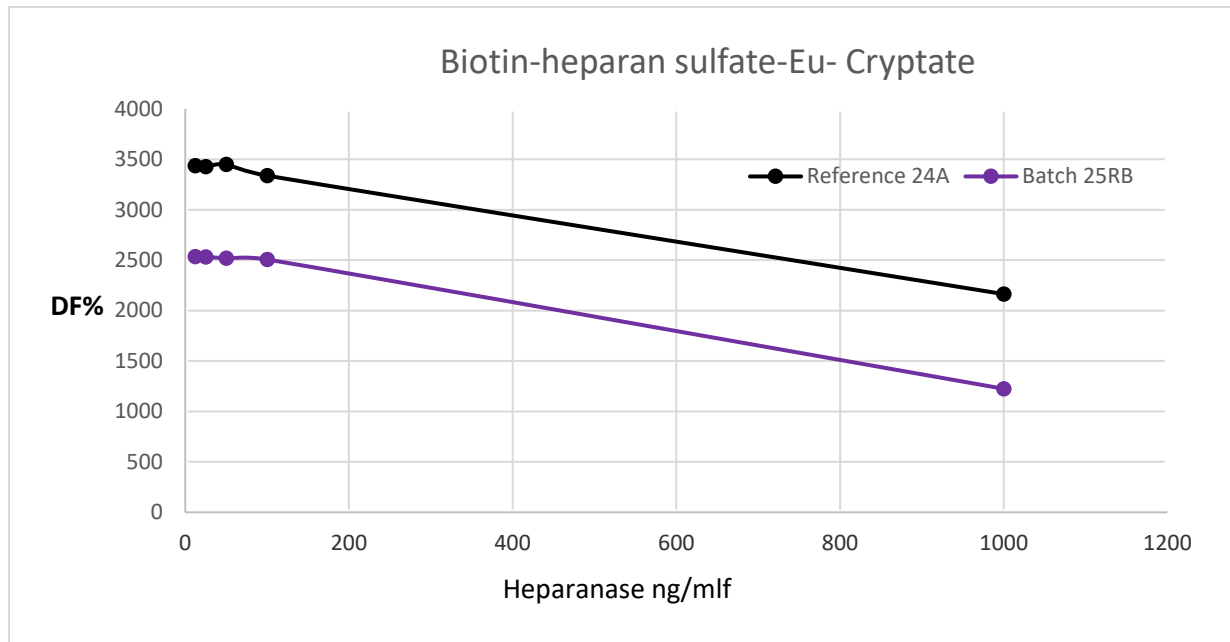
Incubation 30min 37°C

- 10 µL Streptavidin-d2

Incubation 30 min at room temperature. The plate was read on PheraStar Plus under standard conditions.

RESULTS

[Enzyme] ng/mlf		Biot-HS-Cryptate Reference	Biotin-HS-Cryptate Current BATCH	Norm
-	Delta F max%	3 575%	2 499%	>1500%
12.5	DF/DF max%	96%	101%	-
25		96%	101%	-
50		96%	101%	-
100		95%	100%	-
1 000		61%	49%	-
Window		1.59	2.07	-



Biotin-Heparan-sulfate-Eu
Cryptate batch # 25RB

Accepted

Rejected

Operator's initials: MB

Date: 31/07/2018

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.

