

HTRF® Research Reagents

**HTRF CLIP-Red - 20 nmoles**

Product number: SCLPREDE Lot number: 08A

**Product characteristics:**

- Conditioning buffer -
- Appearance Desiccated form
- Storage conditions < -16 °C
- Expiration date 2026.11.14

**Weight indications:**

	Quantity per vial	MW (Da)
Active Moiety (Pro)	20.00 nmoles	N/A

**Please note**

The expiration date of batch 08A has been extended to 2026-11-14 after validation by QC.

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.