

HTRF® Research Reagents

HTRF Biot.Heparan sulfate-Eu - 20K pts
Product number: 61BHSKAB

Lot number: 26A

Product characteristics:

- | | |
|---|---------------------------|
| • Packaging Unit | 5 000 tests/vial |
| • Conditioning buffer | PO4100mM PH7 + 0.1% tween |
| • Appearance | Frozen form |
| • Storage conditions | <-16° °C |
| • Expiration date | 2034.04.22 |
| • Volume per vial | 88.00 µL |
| • Active Moiety Concentration (HS) | 232.70 µg/mL |

Weight indications:

| | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (HS) | 81.00 µg | 11 000 |

Recommended quantity per well*:

| | in 384 well low volume plate (20µL) |
|----------------------|-------------------------------------|
| Active Moiety (HS) | 4.05 ng |
| Number of tests/vial | 5 000 |

*Actual amount per well should be optimized according to assay conditions

Please note

KF can play an essential role in the lanthanide cryptate protection by preventing the action of possible quenchers contained in the assay. KF is mandatory for HTRF assays using Europium cryptate. KF is generally used at a final concentration of 100 to 400 mM, and is added to the conjugate working solutions or dispensed in a separate step, just before the readout.

Once thawed, to avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials (aliquots should not be under 10µL) for storage at -20°C or below.