

**SA-K**

Streptavidin-Eu cryptate - 1 mg - CUSTOM LABELING

Product reference :

**610SAKAH**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 28A  
- Conditioning buffer\_\_\_\_\_ Phosphate 0.1M pH7 + 0.5% BSA +  
0.1% tween  
- Appearance\_\_\_\_\_ Frozen form  
- Storage conditions\_\_\_\_\_ <-16 °C  
- Expiration date\_\_\_\_\_ 2026.01.29  
- Volume per vial\_\_\_\_\_ 4 mL  
- Active Moiety Concentration (SA)\_\_\_\_\_ 250 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (SA)	1.00 mg	60 000

**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.

***Further information as well as a specific application note regarding conjugate-related calculations are available on our HTRF web site: [www.htrf.com/products](http://www.htrf.com/products)***