

**Eu KTBPsuberate**  
Product reference :

Europium cryptate TBP NHS - 0.1 mg  
**65EUSABA**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 25A  
- Conditioning buffer \_\_\_\_\_ Desiccated  
- Appearance \_\_\_\_\_ Desiccated form  
- Storage conditions \_\_\_\_\_ <-16 °C  
- Expiration date \_\_\_\_\_ 2025.04.30

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (K)	0.10 mg	1 449

**PLEASE NOTE**

Rz (305/280) = 3.1

Molar extinction coefficient (305nm) = 30 000 L.Mole<sup>-1</sup>

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