

**KTBP5COOHMP**  
Product reference :

Europium cryptate TBP 5COOH MP - 1 mg  
**65T5MABB**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 07A  
- Conditioning buffer \_\_\_\_\_ Desiccated  
- Appearance \_\_\_\_\_ Desiccated form  
- Storage conditions \_\_\_\_\_ <-16 °C  
- Expiration date \_\_\_\_\_ 2022.09.12

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (K)	1.00 mg	1 523

**PLEASE NOTE**

Rz (320.5/280) = 2.82

Molar extinction coefficient (317 nm) = 31 000 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)***