

## Batch information

**Eu KTBPsuberate**                      Europium cryptate TBP NHS - 1 mg  
Product reference :                      **65EUSABB**

### 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)	1.00 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)***

