

Eu KTBPsuberate
Product reference :

Europium cryptate TBP NHS - 1 mg
65EUSABB

Product characteristics :

- Lot Number _____ 22B
- Conditioning buffer _____ Desiccated
- Appearance _____ Desiccated form
- Storage conditions _____ <-16 °C
- Expiration date _____ 2022.01.10

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (K)	1.00 mg	1 449

PLEASE NOTE

Rz (305/280) = 2.96

Molar extinction coefficient (305nm) = 30 000 L.Mole⁻¹

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com