

Batch information

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

Product characteristics :

- Lot Number	26A
- Conditioning buffer	Desiccated
- Appearance	Desiccated form
- Storage conditions	<-16° °C
- Expiration date	2026.08.16

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⁻¹
Reconstitute with DMSO

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

