

Batch information

STD

HTRF Cereblon Bind. kit - Std

Product reference :

64BDCRBNCDA

Product characteristics :

- Lot Number	01A
- Conditioning buffer	tampon hepes 50mM pH7 + 01% BSA + stabilizer
- Appearance	Frozen form
- Storage conditions	<-60 °C
- Expiration date	2023.06.05

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (LIG)	240.00 nmoles	N/A

PLEASE NOTE

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

