

## Batch information

**STD** Cereblon binding kit - Standard

Product reference : **64BDCRBNCDA**

### Product characteristics :

- Lot Number	02A
- Conditioning buffer	tampon hepes 50mM pH7 + 01% BSA + stabilizer
- Appearance	Frozen form
- Storage conditions	<-60 °C
- Expiration date	2023.11.20

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (LIG)	280.00 µmoles	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](http://www.cisbio.com)*

