

## Batch information

**STD** CD47/SIRP alpha binding kit - Standard

Product reference : **64SIRPCDA**

### Product characteristics :

- Lot Number	02A
- Conditioning buffer	Tampon Hepes 50mM pH 7 + BSA 0.1% + stabilizer
- Appearance	Frozen form
- Storage conditions	< -16 °C
- Expiration date	2024.03.04
- Volume per vial	50.00 µL

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ac)	0.13 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](http://www.cisbio.com)*

