

Batch information

STD

CD47/SIRP alpha binding kit - Standard

Product reference :

64SIRPCDA

Product characteristics :

- Lot Number	03A
- Conditioning buffer	Tampon Hepes 50mM pH 7 + BSA 0.1% + stabilizer
- Appearance	Frozen form
- Storage conditions	< -16 °C
- Expiration date	2024.08.16
- 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

