

STD

CD47/SIRP alpha binding kit - Standard

Product reference :

64SIRPCDA

Product characteristics :

- Lot Number_____ 01A
- Conditioning buffer_____ Tampon Hepes 50mM pH 7 +
BSA 0.1% + stabilizer
- Appearance_____ Frozen form
- Storage conditions_____ < -16 °C
- Expiration date_____ 2023.01.27
- Volume per vial_____ 50 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ac)	0.13 nmoles	N/A

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