

STD

Human STING binding kit - Standard
64BDSTGCDA

Product reference :

Product characteristics :

- Lot Number_____ 02A
- Conditioning buffer_____ Tampon hépés + stabilizer
- Appearance_____ Frozen form
- Storage conditions_____ -60 °C
- Expiration date_____ 2022.05.28
- Volume per vial_____ 50 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	10.00 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

