

STD

Active GLP1 kit calibrator

Product reference :

62GLPCDA

Product characteristics :

- Lot Number_____ 06A
- Conditioning buffer_____ Diluent #4
- Appearance_____ Frozen form
- Storage conditions_____ -60 °C
- Expiration date_____ 2022.07.24
- Volume per vial_____ 50 µL
- Active Moiety Concentration (pep)_____ 56 ng/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (pep)	2.80 ng	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