

Control lysate

YAP total kit control lysate

Product reference :

64YATTDA

Product characteristics :

- Lot Number_____ 09A
- Conditioning buffer_____ Cellular kinase lysis buffer
- Appearance_____ Frozen form
- Storage conditions_____ -60 °C
- Expiration date_____ 2022.11.25
- Volume per vial_____ 150 µL
- Active Moiety Concentration (Ab)_____ 1 X

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	150.00 µl	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

