

Batch information

Control lysate

HTRF FoxO1 p-S256 kit - Ctrl lysate

Product reference :

64FOXTDA

Product characteristics :

- Lot Number	02A
- Conditioning buffer	Lysis Buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.12.03
- Volume per vial	150.00 µL

Weight indications :

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

PLEASE NOTE

The expiration date of batch 02A has been extended to 2024-12-03 after validation by QC.

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

