

Batch information

Alpha-tubulin Standard

Alpha-Tubulin Housekeeping kit - Standard

Product reference :

64ATUBCDA

Product characteristics :

- Lot Number	01A
- Conditioning buffer	N/A
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2023.11.20
- Volume per vial	100.00 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (atu)	3.00 µg	N/A

PLEASE NOTE

The expiration date of batch 01A has been extended to 2023-11-20 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

