

Batch information

Tag-lite labeling medium 5X

SNAP/CLIP lab medium 5X - 100 mL

Product reference :

LABMED

Product characteristics :

- Lot Number	13A
- Conditioning buffer	Stock solution 5 X
- Appearance	Liquid form
- Storage conditions	2-8 ° °C
- Expiration date	2023.01.07
- Volume per vial	100.00 mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	100.00 mL	N/A

PLEASE NOTE

Tag-lite labeling medium 5x must be diluted 1/5 with distilled water under sterile conditions before use

The expiration date og batch 13A has been extended to 2023-01-07 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

