

Batch information

SAH d2 conjugate - Medium

HTRF Methyltransferase kit - 10K pts d2

Product reference :

62SAHDLA

Product characteristics :

- Lot Number	12A
- Conditioning buffer	H2O, 0.1% BSA
- Appearance	Liquid form
- Storage conditions	< -60 °C
- Expiration date	2024.06.23
- Volume per vial	5.50 mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (SAH)	5.50 mL	N/A

PLEASE NOTE

See package insert of #62SAHPEH for use.

The expiration date of batch 12A has been extended to 2024-06-23 after validation by CQ.

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

