

SAH d2 conjugate - Medium

Methyltransferase assay - 10,000 tests d2 conj

Product reference :

62SAHDLA

Product characteristics :

- Lot Number _____ 12A
- Conditioning buffer _____ H2O, 0.1% BSA
- Appearance _____ Liquid form
- Storage conditions _____ < -60 °C
- Expiration date _____ 2022.06.23
- Volume per vial _____ 5.5 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.

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