

Batch information

SAH Standard

Methyltransferase assay - Standard

Product reference :

62SAHCLB

Product characteristics :

- Lot Number	10A
- Conditioning buffer	H2O
- Appearance	Frozen form
- Storage conditions	< -60 °C
- Expiration date	2024.01.11
- Volume per vial	400.00 µL
- Active Moiety Concentration (SAH)	500.00 µM

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (SAH)	400.00 µl	384

PLEASE NOTE

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

