

SAH Standard

Methyltransferase assay standard

Product reference :

62SAHCLB

Product characteristics :

- Lot Number_____ 09A
- Conditioning buffer_____ H2O
- Appearance_____ Frozen form
- Storage conditions_____ < -60 °C
- Expiration date_____ 2022.11.13
- Volume per vial_____ 400 µL
- Active Moiety Concentration (SAH)_____ 500 µM

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (SAH)	0.20 µmoles	384

PLEASE NOTE

The expiration date of batch 09A has been extended to 2022-11-13 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