

## Batch information

**Ultrapure SAM cofactor (2.4 µmol)** Ultrapure SAM cofactor - 2.4 µmol

Product reference : **62SAHZLC**

### Product characteristics :

- Lot Number	15A
- Conditioning buffer	H2O
- Appearance	Frozen form
- Storage conditions	< -60 °C
- Expiration date	2023.11.11
- Volume per vial	1.20 mL
- Active Moiety Concentration (SAM)	2.00 mM

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (SAM)	2.40 µmoles	399

### PLEASE NOTE

The expiration date of batch 15A has been extended to 2023-11-11 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](http://www.cisbio.com)***

