

Ultrapure SAM cofactor (2.4 µmol) Ultrapure SAM cofactor (2.4 µmol)
Product reference : **62SAHZLC**

Product characteristics :

- Lot Number _____ 16A
- Conditioning buffer _____ H₂O
- Appearance _____ Frozen form
- Storage conditions _____ < -60 °C
- Expiration date _____ 2023.02.23
- Volume per vial _____ 1.2 mL
- Active Moiety Concentration (SAM) _____ 2 mM

Weight indications :

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

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