

Ultrapure SAM cofactor (22 µmol) Ultrapure SAM cofactor (22 µmol)
Product reference : **62SAHZLD**

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

PLEASE NOTE

10 vials of 2.4 µmoles (62SAHZLC)

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