

FSK

Forskolin - Adenylyl cyclase activator - 40 µl

Product reference :

62AMYADA

Product characteristics :

- Lot Number_____ 02A
- Conditioning buffer_____ DMSO
- Appearance_____ Liquid form
- Storage conditions_____ 4 °C
- Expiration date_____ 2023.02.15
- Active Moiety Concentration (FSK)_____ 10 mM

Weight indications :

| | Quantity per vial | MW (Da) |
|---------------------|-------------------|---------|
| Active Moiety (FSK) | 40.00 µl | 411 |

PLEASE NOTE

The expiration date of batch 02A has been extended to 2023-02-15 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

