

IBMX

IBMX - Phosphodiesterase inhibitor - 500 µl

Product reference :

62AMXADC

Product characteristics :

- Lot Number_____ 04A
- Conditioning buffer_____ DMSO
- Appearance_____ Liquid form
- Storage conditions_____ 4 °C
- Expiration date_____ 2023.03.19
- Active Moiety Concentration (0)_____ 500 mM

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (0)	500.00 µl	222

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